

PRODUCT INFORMATION



FDA-Approved Drugs Screening Library

Item No. 23538 • Batch No. 0646892

Panels are routinely re-evaluated to include new catalog introductions as the research evolves.

Page 1 of 32

Plate	Well	Contents	Item Number	Primary Target*
1	A1	Unused		
1	A2	Acetaminophen	10024	
1	A3	Treprostinil	10162	DP ₁ , EP ₂ , IP Receptors
1	A4	PKC 412	10459	Multiple Kinases
1	A5	Paclitaxel	10461	Tubulin
1	A6	Erlotinib	10483	EGFR
1	A7	Capecitabine	10487	Thymidylate Synthase
1	A8	Bupropion (hydrochloride)	10488	DAT, NET
1	A9	Candesartan Cilexetil	10489	AT ₁ Receptors
1	A10	Atorvastatin (calcium salt hydrate)	10493	HMG-CoA Reductase
1	A11	PLX4032	10618	B-Raf, B-Raf ^{V600E}
1	A12	Unused		
1	B1	Unused		
1	B2	Olaparib	10621	PARP1, PARP2
1	B3	Guanabenz (hydrochloride)	10851	α ₂ -Adrenergic Receptors
1	B4	Dexamethasone	11015	Glucocorticoid Receptors
1	B5	all-trans Retinoic Acid	11017	RAR
1	B6	BIBF 1120	11022	FGFR, PDGFR, VEGFR
1	B7	6-ECDCA	11031	FXR
1	B8	Nifedipine	11106	Ca _v 1 Channels
1	B9	Diphenhydramine (hydrochloride)	11158	Histamine H ₁ , Muscarinic M ₁₋₅ Receptors
1	B10	5-Azacytidine	11164	DNA Methyltransferase
1	B11	Decitabine	11166	DNA Methyltransferase
1	B12	Unused		
1	C1	Unused		
1	C2	Sodium 4-Phenylbutyrate	11323	Aberrant Protein Folding
1	C3	Actinomycin D	11421	DNA (Intercalating Agent)
1	C4	Lapatinib	11493	EGFR, HER2
1	C5	Ponatinib	11494	Bcr-Abl
1	C6	Dasatinib	11498	Abl, LCK, Src
1	C7	Venlafaxine (hydrochloride)	11567	DAT, NET, SERT
1	C8	Letrozole	11568	Aromatase
1	C9	Bexarotene	11571	RXR
1	C10	Dapagliflozin	11574	SGLT
1	C11	Canagliflozin	11575	SGLT
1	C12	Unused		

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/04/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 2 of 32

Plate	Well	Contents	Item Number	Primary Target*
1	D1	<i>Unused</i>		
1	D2	Eprosartan (mesylate)	11578	AT ₁ Receptors
1	D3	Glipizide	11579	K _{ATP} Channels
1	D4	Temsirolimus	11590	FKBP-12-mTOR Interaction
1	D5	Fluconazole	11594	Ergosterol Biosynthesis
1	D6	MDV 3100	11596	Androgen Receptors
1	D7	Everolimus	11597	FKBP-12-mTOR Interaction
1	D8	Tofacitinib (citrate)	11598	JAK1, JAK2, JAK3
1	D9	Olmesartan Medoxomil	11614	AT ₁ Receptors
1	D10	Telmisartan	11615	AT ₁ Receptors
1	D11	Tamoxifen (citrate)	11629	Estrogen Receptors
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	Natamycin	11634	Ergosterol Biosynthesis
1	E3	5-Fluorocytosine	11635	DNA Synthesis
1	E4	Amphotericin B	11636	Ergosterol Biosynthesis
1	E5	Docetaxel (hydrate)	11637	Tubulin
1	E6	Brinzolamide	11660	CAII, CAIV
1	E7	Gemcitabine	11690	DNA (Intercalating Agent)
1	E8	Bosentan	11731	ET _A Receptors
1	E9	Atazanavir	11733	HIV-1 Protease
1	E10	Vinblastine (sulfate)	11762	Tubulin
1	E11	Vincristine (sulfate)	11764	Tubulin
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	Vitamin D ₂	11791	Calcium Levels
1	F3	Oxytocin (acetate)	11799	OXTR
1	F4	Artemether	11815	
1	F5	Rosiglitazone (maleate)	11884	PPAR γ
1	F6	Acarbose	11885	α -Glucosidase
1	F7	Olanzapine	11937	Various CNS Receptors
1	F8	Irbesartan	11952	AT ₁ Receptors
1	F9	Fingolimod	11975	S1P
1	F10	(S)-Pramipexole (hydrochloride)	11981	Dopamine Receptors
1	F11	Anastrozole	11987	Aromatase
1	F12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 3 of 32

Plate	Well	Contents	Item Number	Primary Target*
1	G1	<i>Unused</i>		
1	G2	Olopatadine (hydrochloride)	11999	Histamine H ₁ Receptors
1	G3	Atropine	12008	Muscarinic Receptors
1	G4	Haloperidol	12014	Dopamine Receptors
1	G5	Bosutinib	12030	Abl, EPHB2, Src
1	G6	Clozapine	12059	5-HT, Histamine, α ₁ -Adrenergic Receptors
1	G7	1-hydroxy Vitamin D ₂	12077	Vitamin D Receptor
1	G8	Cladribine	12085	Lymphocyte DNA Synthesis
1	G9	(R)-Crizotinib	12087	ALK, c-MET
1	G10	Cyclosporin A	12088	Cyclophilin D
1	G11	Glimepiride	12090	K _{ATP} Channels
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	Epirubicin (hydrochloride)	12091	DNA (Intercalating Agent)
1	H3	Etoposide	12092	Topoisomerase II
1	H4	Pazopanib	12097	Multiple Kinases
1	H5	Prostaglandin E ₁	13010	EP ₁ Receptors
1	H6	Valproic Acid (sodium salt)	13033	HDAC
1	H7	Streptozotocin	13104	DNA (Alkylating Agent)
1	H8	Oxaliplatin	13106	DNA (Cross-linking Agent)
1	H9	Pimecrolimus	13107	Calcineurin, Prolyl Isomerase
1	H10	Carboplatin	13112	DNA (Cross-linking Agent)
1	H11	Cidofovir	13113	DNA Polymerase
1	H12	<i>Unused</i>		
2	A1	<i>Unused</i>		
2	A2	Metformin (hydrochloride)	13118	Multiple Targets
2	A3	Cisplatin	13119	DNA (Cross-linking Agent)
2	A4	Imatinib (mesylate)	13139	Multiple Kinases
2	A5	Sunitinib (malate)	13159	FLK1, FLT3, PDGFRβ
2	A6	Donepezil	13245	AChE
2	A7	Eltrombopag	13247	Thrombopoietin Receptors
2	A8	(-)-Sitagliptin (phosphate)	13252	DPP-4
2	A9	Panobinostat	13280	HDAC
2	A10	Itraconazole	13288	Ergosterol biosynthesis
2	A11	MCI-186	13320	Oxidative Stress
2	A12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 4 of 32

Plate	Well	Contents	Item Number	Primary Target*
2	B1	<i>Unused</i>		
2	B2	Rapamycin	13346	mTORC1
2	B3	Milrinone	13357	PDE3
2	B4	Piroxicam	13368	COX, NSAID
2	B5	Trandolapril	13522	ACE
2	B6	(±)-Warfarin	13566	Vitamin K
2	B7	Topiramate	13623	Multiple Targets
2	B8	Lansoprazole	13627	H ⁺ /K ⁺ ATPase
2	B9	Risperidone	13629	Multiple Targets
2	B10	Alendronate (sodium hydrate)	13642	Osteoclasts
2	B11	Alfuzosin (hydrochloride)	13648	α ₁ -Adrenergic Receptors
2	B12	<i>Unused</i>		
2	C1	<i>Unused</i>		
2	C2	Adapalene	13655	RAR
2	C3	(±)-Clopidogrel (hydrochloride)	13657	P2Y ₁₂ Receptors
2	C4	Nitazoxanide	13692	Pyruvate:Ferredoxin Oxidoreductase
2	C5	Axitinib	13813	VEGFR
2	C6	Misoprostol	13820	EP ₂ , EP ₃ Receptors
2	C7	Entecavir (hydrate)	13831	HBV Reverse Transcriptase
2	C8	Cyclophosphamide (hydrate)	13849	DNA (Alkylating Agent)
2	C9	Ganciclovir	13853	DNA Polymerase
2	C10	Lopinavir	13854	HIV-1 Protease
2	C11	Ritonavir	13872	HIV-1 Protease
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	Tenofovir	13874	HIV-1 Reverse Transcriptase
2	D3	Valganciclovir (hydrochloride)	13875	DNA Polymerase
2	D4	Bleomycin (sulfate)	13877	
2	D5	Dextromethorphan (hydrobromide hydrate)	13950	σ-Opioid Receptors
2	D6	Methotrexate (hydrate)	13960	Folate Metabolism
2	D7	Timolol (maleate)	13974	β-Adrenergic Receptors
2	D8	Pirfenidone	13986	
2	D9	Azathioprine	13987	
2	D10	Mycophenolate Mofetil	13988	IMPDH
2	D11	Sildenafil Citrate	14008	PDE5
2	D12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 5 of 32

Plate	Well	Contents	Item Number	Primary Target*
2	E1	<i>Unused</i>		
2	E2	Prostaglandin E ₂	14010	EP Receptors
2	E3	Tadalafil	14024	PDE5
2	E4	Ingenol	14031	PKC
2	E5	Propylthiouracil	14069	Thyroid Peroxidase
2	E6	Scopolamine (hydrobromide)	14108	Muscarinic Receptors
2	E7	L-Thyroxine	14116	Thyroid Receptors
2	E8	Caffeine	14118	Multiple Targets
2	E9	Clofarabine	14125	Deoxycytidine Kinase, DNA Polymerase
2	E10	Febuxostat	14127	Xanthine Oxidase
2	E11	Topotecan (hydrochloride)	14129	Topoisomerase I
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	Quetiapine (hemifumarate)	14152	
2	F3	Entacapone	14153	Catechol O-methyltransferase
2	F4	Floxuridine	14154	Multiple Targets
2	F5	Polymyxin B (sulfate)	14157	LPS
2	F6	Daunorubicin (hydrochloride)	14159	Topoisomerase
2	F7	Acyclovir	14160	DNA Polymerase
2	F8	Temozolomide	14163	DNA (Alkylating Agent)
2	F9	Idarubicin (hydrochloride)	14176	Topoisomerase II
2	F10	Valsartan	14178	AT ₂ Receptors
2	F11	Irinotecan (hydrochloride hydrate)	14180	Topoisomerase I
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	Memantine (hydrochloride)	14184	NMDA Receptors
2	G3	Chloroquine (phosphate)	14194	Autophagosome-Lysosome Fusion
2	G4	Bicalutamide	14250	Androgen Receptors
2	G5	Fludarabine Phosphate	14251	
2	G6	Flumazenil	14252	GABA _A Receptors
2	G7	<i>Unused</i>		
2	G8	Rivastigmine (tartrate)	14270	AChE, BChE
2	G9	Prasugrel	14278	P2Y ₁₂ Receptors
2	G10	Ciprofloxacin (hydrochloride hydrate)	14286	Topoisomerase II, Topoisomerase IV
2	G11	<i>Unused</i>		
2	G12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 6 of 32

Plate	Well	Contents	Item Number	Primary Target*
2	H1	<i>Unused</i>		
2	H2	Verapamil (hydrochloride)	14288	L-type Calcium Channels
2	H3	Gatifloxacin	14290	DNA Gyrase
2	H4	(S)-Duloxetine (hydrochloride)	14317	NET, SERT
2	H5	Dantrolene (sodium salt)	14326	Calcium Release
2	H6	Tetracycline (hydrochloride)	14328	
2	H7	<i>Unused</i>		
2	H8	Thioridazine (hydrochloride)	14400	Multiple Targets
2	H9	A-771726	14404	DHODH
2	H10	Amiloride (hydrochloride) (hydrate)	14409	Potassium
2	H11	Efavirenz	14412	Reverse Transcriptase
2	H12	<i>Unused</i>		
3	A1	<i>Unused</i>		
3	A2	5-Fluorouracil	14416	Thymidylate Synthase
3	A3	Ampicillin (sodium salt)	14417	Transpeptidase
3	A4	Fluoxetine (hydrochloride)	14418	SERT
3	A5	Doxycycline (hydrate)	14422	Bacterial 30S Ribosomal Subunit
3	A6	Rifampicin	14423	RNA Polymerase
3	A7	Darifenacin (hydrobromide)	14424	Muscarinic M ₁₋₅ Receptors
3	A8	Teniposide	14425	Topoisomerase II
3	A9	Minocycline (hydrochloride hydrate)	14454	Bacterial 30S Ribosomal Subunit
3	A10	Carbamoylcholine (chloride)	14486	Muscarinic Receptors, nAChRs
3	A11	(+)-Pilocarpine (hydrochloride)	14487	Muscarinic M ₁₋₃ Receptors
3	A12	<i>Unused</i>		
3	B1	<i>Unused</i>		
3	B2	Citalopram (hydrobromide)	14572	SERT
3	B3	Nimodipine	14573	Ca _v 1.2 Channels
3	B4	Diazoxide	14576	SUR1/K _{IR} 6.2
3	B5	Riluzole (hydrochloride)	14577	Glutamate Neurotransmission
3	B6	9- <i>cis</i> -Retinoic Acid	14587	RAR, RXR
3	B7	Deferoxamine (mesylate)	14595	Iron
3	B8	Tobramycin	14596	Bacterial 30S Ribosomal Subunit, Bacterial 50S Ribosomal Subunit
3	B9	Bromocriptine (mesylate)	14598	Various CNS Receptors
3	B10	Sumatriptan (succinate)	14600	5-HT _{1B} , 5-HT _{1D} , 5-HT _{1F} Receptors
3	B11	Mecamylamine (hydrochloride)	14602	nAChRs (α3β4 Subunit)
3	B12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 7 of 32

Plate	Well	Contents	Item Number	Primary Target*
3	C1	<i>Unused</i>		
3	C2	Meclizine (hydrochloride)	14615	Histamine H ₁ Receptors
3	C3	Dorzolamide (hydrochloride)	14616	Carbonic Anhydrase
3	C4	Bumetanide	14630	NKCC1
3	C5	Homoharringtonine	14631	Human 80S Ribosome
3	C6	Dexrazoxane	14632	Iron, Topoisomerase 2A
3	C7	Clemastine (fumarate)	14637	Histamine H ₁ , Muscarinic Receptors
3	C8	Maraviroc	14641	CCR5
3	C9	Lenalidomide	14643	CRBN
3	C10	Vandetanib	14706	Multiple Kinases
3	C11	Dimethyl fumarate	14714	Nrf2
3	C12	<i>Unused</i>		
3	D1	<i>Unused</i>		
3	D2	Posaconazole	14737	CYP51
3	D3	Abacavir (sulfate)	14746	Reverse Transcriptase
3	D4	Ceftazidime (hydrate)	14828	PBP3
3	D5	Moxifloxacin (hydrochloride)	14830	DNA Gyrase
3	D6	Gemfibrozil	14835	PPAR α , PPAR γ
3	D7	Amlodipine	14838	Ca _v Channels
3	D8	Sertraline (hydrochloride)	14839	SERT
3	D9	Mitoxantrone (hydrochloride)	14842	DNA (Intercalating Agent), Topoisomerase II, HIV Integrase
3	D10	Busulfan	14843	DNA (Alkylating Agent)
3	D11	Leflunomide	14860	AhR, DHODH
3	D12	<i>Unused</i>		
3	E1	<i>Unused</i>		
3	E2	Sildenafil	14866	α_{1A} -Adrenergic Receptors
3	E3	Aprepitant	14867	NK1
3	E4	Loperamide (hydrochloride)	14875	μ -Opioid Receptors
3	E5	Pentostatin	14878	ADA
3	E6	Omeprazole	14880	H ⁺ /K ⁺ ATPase
3	E7	Meloxicam	14906	COX-2
3	E8	Rasagiline	14917	MAO-B
3	E9	Vardenafil (hydrochloride hydrate)	14930	PDE5, PDE6
3	E10	Finasteride	14938	5 α -Reductase Type II
3	E11	Rabeprazole (sodium salt)	14939	H ⁺ /K ⁺ ATPase
3	E12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 8 of 32

Plate	Well	Contents	Item Number	Primary Target*
3	F1	<i>Unused</i>		
3	F2	Imiquimod	14956	TLR7
3	F3	Stavudine	14975	Reverse Transcriptase
3	F4	Probenecid	14981	OATs, MRP1, TRPV2
3	F5	<i>Unused</i>		
3	F6	Metyrapone	14994	CYP450
3	F7	Paroxetine (hydrochloride)	14998	SERT
3	F8	Ondansetron (hydrochloride hydrate)	14999	5-HT ₃ Receptors
3	F9	Azithromycin	15004	Bacterial 50S Ribosomal Subunit
3	F10	Bazedoxifene (acetate)	15005	ER α , ER β
3	F11	Clindamycin	15006	Bacterial 50S Ribosomal Subunit
3	F12	<i>Unused</i>		
3	G1	<i>Unused</i>		
3	G2	Doxorubicin (hydrochloride)	15007	DNA (Intercalating Agent), Topoisomerase II
3	G3	Exemestane	15008	Steroidal Aromatase
3	G4	Glyburide (potassium salt)	15009	SUR1/K _{IR} 6.2
3	G5	(\pm)-Ibutilide (hemifumarate)	15011	I _{Kr}
3	G6	Linezolid	15012	Bacterial 50S Ribosomal Subunit
3	G7	Methylprednisolone	15013	Glucocorticoid, Mineralocorticoid Receptors
3	G8	Miglitol	15014	α -Glucosidase
3	G9	Prazosin (hydrochloride)	15023	α_1 -Adrenergic, MT ₃ Receptors
3	G10	Sulfasalazine	15025	
3	G11	Tigecycline	15026	Bacterial 30S Ribosomal Subunit
3	G12	<i>Unused</i>		
3	H1	<i>Unused</i>		
3	H2	Tolterodine (tartrate)	15027	Muscarinic M ₁₋₅ Receptors
3	H3	Varenicline (tartrate)	15030	nAChRs (α 7)
3	H4	Ziprasidone (hydrochloride hydrate)	15031	Various CNS Receptors
3	H5	Cilostazol	15035	PDE3A
3	H6	Dofetilide	15045	I _{Kr}
3	H7	R-(-)-Deprenyl (hydrochloride)	15046	MAO-B
3	H8	Trifluoperazine (hydrochloride)	15068	α_{1A} -Adrenergic, Dopamine D ₂ Receptors
3	H9	Nevirapine	15117	Reverse Transcriptase
3	H10	Zanamivir	15123	Neuraminidase
3	H11	Roflumilast	15141	PDE4B, PDE4D
3	H12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 9 of 32

Plate	Well	Contents	Item Number	Primary Target*
4	A1	<i>Unused</i>		
4	A2	Nelfinavir (mesylate)	15144	HIV-1 Protease
4	A3	Ivacaftor	15145	CFTR
4	A4	Abiraterone Acetate	15148	CYP17A1
4	A5	Indinavir (sulfate)	15150	HIV-1 Protease
4	A6	Ketoconazole	15212	Sterol 14 α Demethylase
4	A7	Amiodarone (hydrochloride)	15213	hERG
4	A8	Clotrimazole	15278	Sterol 14 α Demethylase
4	A9	CAL-101	15279	PI3K p110 δ
4	A10	Minoxidil	15302	COX-1
4	A11	Disulfiram	15303	Aldehyde Dehydrogenase
4	A12	<i>Unused</i>		
4	B1	<i>Unused</i>		
4	B2	Captopril	15313	ACE
4	B3	Auranofin	15316	Thioredoxin Reductase
4	B5	Vancomycin (hydrochloride)	15327	Transpeptidation
4	B6	Amprenavir	15369	HIV Protease
4	B7	PSI-7977	15402	DNA Replication
4	B8	Amikacin	15405	Bacterial 30S Ribosomal Subunit
4	B9	Budesonide	15407	Glucocorticoid Receptors
4	B10	Pitavastatin (calcium salt)	15414	HMG-CoA Reductase
4	B11	Carvedilol	15418	β -Adrenergic Receptors
4	B12	<i>Unused</i>		
4	C1	<i>Unused</i>		
4	C2	Miconazole (nitrate)	15420	Sterol 14 α Demethylase
4	C3	Ticagrelor	15425	P2Y ₁₂ Receptors
4	C4	Brimonidine (tartrate)	15426	α_2 -Adrenergic Receptors
4	C5	Apixaban	15427	Factor Xa
4	C6	Lamotrigine	15428	Na _v Channels
4	C7	Oxaprozin	15476	COX-1, COX-2
4	C8	Zidovudine	15492	Reverse Transcriptase
4	C9	Fidaxomicin	15503	RNA Polymerase
4	C10	Naltrexone (hydrochloride)	15520	Opioid Receptors
4	C11	Paliperidone	15556	Dopamine D ₂ , D ₃ Receptors
4	C12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 10 of 32

Plate	Well	Contents	Item Number	Primary Target*
4	D1	<i>Unused</i>		
4	D2	Ramipril	15558	ACE
4	D3	Dexmedetomidine (hydrochloride)	15581	α_2 -Adrenergic Receptors
4	D4	Dobutamine (hydrochloride)	15582	β -Adrenergic Receptors
4	D5	Formoterol (hemifumarate hydrate)	15584	β_2 -Adrenergic Receptors
4	D6	Isoproterenol (hydrochloride)	15592	β -Adrenergic Receptors
4	D7	Naloxone (hydrochloride)	15594	Opioid Receptors
4	D8	Vecuronium (bromide)	15603	nAChRs
4	D9	Ranolazine (hydrochloride)	15604	I_{Na}
4	D10	Daptomycin	15615	Bacterial Cell Membrane
4	D11	Eplerenone	15616	Mineralocorticoid Receptors
4	D12	<i>Unused</i>		
4	E1	<i>Unused</i>		
4	E2	Fluvoxamine (maleate)	15617	Various CNS Receptors
4	E3	Voriconazole	15633	Sterol 14 α Demethylase
4	E4	Rucaparib	15643	PARP1
4	E5	Dehydroepiandrosterone	15728	
4	E6	Tiotropium (bromide)	15773	Muscarinic M ₁₋₃ Receptors
4	E7	6-Thioguanine	15774	RNA Polymerase
4	E8	Carmustine	15775	DNA (Alkylating Agent)
4	E9	Cefazolin (sodium salt)	15776	Penicillin-binding Proteins
4	E10	Metaxalone	15777	
4	E11	Darunavir	15866	HIV-1 Protease
4	E12	<i>Unused</i>		
4	F1	<i>Unused</i>		
4	F2	Ivabradine (hydrochloride)	15868	HCN Channel
4	F3	Zolmitriptan	15871	5-HT _{1B} , 5-HT _{1D} Receptors
4	F4	Progesterone	15876	Progesterone Receptors
4	F5	Amitriptyline (hydrochloride)	15881	NET, SERT
4	F6	Clomipramine (hydrochloride)	15884	NET, SERT
4	F7	Cyclobenzaprine (hydrochloride)	15885	
4	F8	Doxepin (hydrochloride)	15888	NET, SERT
4	F9	Imipramine (hydrochloride)	15890	NET, SERT
4	F10	Maprotiline (hydrochloride)	15892	NET
4	F11	Nortriptyline (hydrochloride)	15904	NET, SERT
4	F12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number	Primary Target*
4	G1	<i>Unused</i>		
4	G2	Trimipramine (maleate)	15921	SERT
4	G3	Caspofungin (acetate)	15923	1,3-β-D-Glucan Synthase
4	G4	Edrophonium (chloride)	15928	AChE
4	G5	Clonidine (hydrochloride)	15949	α ₂ -Adrenergic Receptors
4	G6	Dutasteride	15956	5α-Reductase
4	G7	Metolazone	15987	Na ⁺ /Cl ⁺ Cotransporter
4	G8	Lofexidine (hydrochloride)	15989	α ₂ -Adrenergic Receptors
4	G9	Cefoxitin (sodium salt)	15990	Penicillin-binding Proteins
4	G10	Salmeterol (xinafoate)	16009	β ₂ -Adrenergic Receptors
4	G11	Ciclopirox	16021	HIF-PH2
4	G12	<i>Unused</i>		
4	H1	<i>Unused</i>		
4	H2	Cefotaxime (sodium salt)	16040	β-Lactamase
4	H3	Enalapril (maleate)	16041	ACE
4	H4	Cinacalcet (hydrochloride)	16042	CaSR
4	H5	Rivaroxaban	16043	Factor Xa
4	H6	Meropenem (hydrate)	16068	Bacterial Cell Wall Synthesis
4	H7	Cytarabine	16069	DNA Polymerase
4	H8	Oseltamivir (phosphate)	16070	Neuraminidase
4	H9	Raltegravir (potassium salt)	16071	HIV-1 Integrase
4	H10	Cefpodoxime	16075	Bacterial Cell Wall Synthesis
4	H11	Clomiphene (citrate)	16087	Estrogen Receptors
4	H12	<i>Unused</i>		
5	A1	<i>Unused</i>		
5	A2	(-)-Apomorphine (hydrochloride hydrate)	16094	Dopamine Receptors
5	A3	Chlorpromazine (hydrochloride)	16129	Various CNS Receptors
5	A4	Rifaximin	16131	PXR
5	A5	Procarbazine (hydrochloride)	16133	DNA Methylation
5	A6	Phentolamine (mesylate)	16135	α-Adrenergic Receptors
5	A7	Sotalol (hydrochloride)	16136	β-Adrenergic Receptors
5	A8	Phenoxybenzamine (hydrochloride)	16211	α-Adrenergic Receptors
5	A9	Chlorthalidone	16213	Na ⁺ /Cl ⁻ Symporter
5	A10	Benztropine (mesylate)	16214	Muscarinic M ₁ Receptors
5	A11	Pimozide	16222	Dopamine Receptors
5	A12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number	Primary Target*
5	B1	<i>Unused</i>		
5	B2	ABT-199	16233	Bcl-2
5	B3	AZD 9291	16237	EGFR ^{T790M}
5	B4	LDE225 (phosphate)	16263	Smoothened
5	B5	PD 0332991 (hydrochloride)	16273	Cdk4, Cdk6
5	B6	Ibrutinib	16274	BTK
5	B7	Trametinib	16292	MEK1, MEK2
5	B8	Ezetimibe	16331	NPC1L1
5	B9	Rifabutin	16468	Bacterial DNA-dependent RNA Polymerase
5	B10	Danazol	16471	Androgen Receptors
5	B11	Pyrimethamine	16472	DHFR
5	B12	<i>Unused</i>		
5	C1	<i>Unused</i>		
5	C2	Trimethoprim	16473	DHFR
5	C3	Reserpine	16474	VMAT1, VMAT2
5	C4	Rizatriptan (benzoate salt)	16475	5-HT _{1B} , 5-HT _{1D} Receptors
5	C5	Tizanidine (hydrochloride)	16477	α ₂ -Adrenergic Receptors
5	C6	Promethazine (hydrochloride)	16478	Histamine H ₁ Receptors
5	C7	Erythromycin	16486	Bacterial 50S Ribosomal Subunit
5	C8	Glycopyrrolate	16498	Muscarinic Receptors
5	C9	(±)-Nicotine	16535	nAChRs
5	C10	Tropicamide	16606	Muscarinic M ₄ Receptors
5	C11	Bupivacaine	16618	Sodium Channels
5	C12	<i>Unused</i>		
5	D1	<i>Unused</i>		
5	D2	Melphalan	16665	DNA (Alkylating Agent)
5	D3	(-)-Norepinephrine (bitartrate hydrate)	16673	α ₁ -Adrenergic, α ₂ -Adrenergic, β ₁ -Adrenergic Receptors
5	D4	Baricitinib	16707	JAK1, JAK2
5	D5	Palonosetron (hydrochloride)	16752	5-HT ₃ Receptors
5	D6	Deferasirox	16753	Iron
5	D7	Ribavirin	16757	Viral Polymerases
5	D8	Fluprostenol isopropyl ester	16769	FP Receptors
5	D9	IPI-145	16800	PI3Kγ, PI3Kδ
5	D10	Latanoprost	16812	FP Receptors
5	D11	17-phenyl trinor Prostaglandin F _{2α} ethyl amide	16820	FP Receptors
5	D12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 13 of 32

Plate	Well	Contents	Item Number	Primary Target*
5	E1	<i>Unused</i>		
5	E2	Lisinopril	16833	ACE
5	E3	Emtricitabine	16879	HBV, HIV-1 Reverse Transcriptases
5	E4	DL- α -Difluoromethylornithine (hydrochloride hydrate)	16889	Ornithine Decarboxylase
5	E5	Tenofovir Disoproxil (fumarate)	16922	HBV, HIV-1 Reverse Transcriptases
5	E6	Desloratadine	16931	Histamine H ₁ Receptors
5	E7	Doripenem (hydrate)	16934	Bacterial Cell Wall Synthesis
5	E8	Ranitidine (hydrochloride)	16939	Histamine H ₂ Receptors
5	E9	Dabrafenib	16989	Raf
5	E10	Binimetinib	16996	MEK1, MEK2
5	E11	Romidepsin	17130	Class I HDACs
5	E12	<i>Unused</i>		
5	F1	<i>Unused</i>		
5	F2	Dabigatran etexilate	17131	Thrombin
5	F3	ARN-509	17132	Androgen Receptors
5	F4	Angiotensin II (human) (acetate)	17150	AT ₁ , AT ₂ Receptors
5	F5	Moxidectin	17165	GABA _A Receptors
5	F6	Cefixime	17176	Bacterial Cell Wall Synthesis
5	F7	1-Deoxygalactonojirimycin (hydrochloride)	17179	α -Galactosidase A
5	F8	Colistin Methanesulfonate (sodium salt)	17180	Bacterial Cell Wall
5	F9	(\pm)-Atenolol	17250	β_1 -Adrenergic Receptors
5	F10	Betaine Aldehyde (chloride)	17270	Homocysteine
5	F11	Furosemide	17273	NKCC1, NKCC2
5	F12	<i>Unused</i>		
5	G1	<i>Unused</i>		
5	G2	Mirabegron	17319	β_3 -Adrenergic Receptors
5	G3	Solifenacin (succinate)	17320	Muscarinic M ₁₋₃ Receptors
5	G4	Esomeprazole Magnesium	17326	H ⁺ /K ⁺ ATPase
5	G5	Verteporfin	17334	Endothelium
5	G6	Oxcarbazepine	17340	Calcium, Sodium, Potassium Channels
5	G7	Chlorhexidine (hydrochloride)	17343	Bacterial Cell Wall
5	G8	Amoxapine	17347	NET, SERT, Various CNS Receptors
5	G9	Desmopressin (trifluoroacetate salt)	17348	Vasopressin V _{1a} , V _{1b} , V ₂ Receptors
5	G10	Midodrine (hydrochloride)	17349	α_1 -Adrenergic Receptors
5	G11	Empagliflozin	17375	SGLT2
5	G12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 14 of 32

Plate	Well	Contents	Item Number	Primary Target*
5	H1	<i>Unused</i>		
5	H2	Felodipine	17535	Ca _v 1 Channels
5	H3	Isradipine	17536	Ca _v 1 Channels
5	H4	Nicardipine (hydrochloride)	17537	Ca _v 1 Channels
5	H5	Carfilzomib	17554	20S Proteasome β5 Subunit
5	H6	Letermovir	17556	CMV DNA Terminase Complex
5	H7	Galantamine	17559	nAChRs
5	H8	Ifosfamide	17562	DNA (Cross-linking Agent)
5	H9	Enalaprilat (hydrate)	17603	ACE
5	H10	Fenoldopam (mesylate)	17629	Dopamine D _{1A} , D _{1B} Receptors
5	H11	LEE011	17666	Cdk4, Cdk6
5	H12	<i>Unused</i>		
6	A1	<i>Unused</i>		
6	A2	Stiripentol	17781	GABA _A Receptors
6	A3	Atracurium (besylate)	17796	Muscarinic Receptors, nAChRs
6	A4	Elvitegravir	17798	HIV-1 Integrase
6	A5	Cevimeline (hydrochloride)	17802	Muscarinic Receptors
6	A6	Fomepizole	17851	Alcohol Dehydrogenase
6	A7	Articaine (hydrochloride)	17883	Na _v Channels
6	A8	Chlorothiazide	17909	
6	A9	Hydroxychloroquine (sulfate)	17911	
6	A10	Nitisinone	17924	4-Hydroxyphenylpyruvate Dioxygenase
6	A11	Felbamate	18000	
6	A12	<i>Unused</i>		
6	B1	<i>Unused</i>		
6	B2	Micafungin (sodium salt)	18009	1,3-β-D-Glucan Synthase
6	B3	Triamcinolone acetone	18026	Glucocorticoid Receptors
6	B4	Oxytetracycline	18076	Amino-Acyl tRNA
6	B5	Epinastine (hydrochloride)	18136	Histamine H ₁ Receptors
6	B6	Riboflavin 5'-monophosphate (sodium salt hydrate)	18167	Corneal Collagen Cross-Linking
6	B7	Dipyridamole	18189	PDE5A
6	B8	Fexofenadine (hydrochloride)	18191	Histamine H ₁ Receptors
6	B9	Tolmetin (sodium salt hydrate)	18195	NSAID
6	B10	Iloprost	18215	IP, EP ₁ Receptors
6	B11	Prostaglandin I ₂ (sodium salt)	18220	Platelets
6	B12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 15 of 32

Plate	Well	Contents	Item Number	Primary Target*
6	C1	<i>Unused</i>		
6	C2	Primaquine (phosphate)	18223	
6	C3	Ciclesonide	18227	Glucocorticoid Receptors
6	C4	(R)-Lansoprazole	18235	H ⁺ /K ⁺ ATPase
6	C5	MLN9708	18386	20S Proteasome
6	C6	Neratinib	18404	EGFR, HER2
6	C7	Vitamin B ₁₂ (hydrate)	18425	Coenzyme B ₁₂
6	C8	Iopromide	18470	Radiographic Contrast Agent
6	C9	Regorafenib	18498	Multiple Kinases
6	C10	R788 (sodium salt)	18501	Tyrosine Kinases
6	C11	Apremilast	18502	PDE4
6	C12	<i>Unused</i>		
6	D1	<i>Unused</i>		
6	D2	4-Aminopyridine	18511	Potassium Channels
6	D3	Lamivudine	18514	HIV-1 Reverse Transcriptase
6	D4	CH5424802	18516	ALK
6	D5	Ledipasvir	18519	HCV NS5A
6	D6	(±)-Baclofen	18600	GABA Receptors
6	D7	Betaxolol (hydrochloride)	18625	β ₁ -Adrenergic Receptors
6	D8	Doxazosin (mesylate)	18633	α ₁ -Adrenergic Receptors
6	D9	Adefovir	18650	HBV DNA Polymerase
6	D10	Mitotane	18651	
6	D11	Balsalazide	18680	
6	D12	<i>Unused</i>		
6	E1	<i>Unused</i>		
6	E2	Pentoxifylline	18720	Erythrocytes
6	E3	Oxiconazole (nitrate)	18724	Ergosterol Biosynthesis
6	E4	Cimetidine	18743	Histamine H ₂ Receptors
6	E5	Regadenoson	18814	
6	E6	Nizatidine	18861	Histamine H ₂ Receptors
6	E7	Ceftriaxone (sodium salt hydrate)	18866	Bacterial Cell Wall Synthesis
6	E8	Chlorzoxazone	18869	
6	E9	Rufinamide	18870	
6	E10	Mebendazole	18872	Tubulin
6	E11	Droperidol	18884	Dopamine Receptors
6	E12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 16 of 32

Plate	Well	Contents	Item Number	Primary Target*
6	F1	<i>Unused</i>		
6	F2	Methsuximide	19058	Calcium Channels
6	F3	Flavoxate (hydrochloride)	19096	Ca _v 1.2 Channels, Muscarinic Receptors
6	F4	Diflunisal	19187	NSAID
6	F5	Amoxicillin (hydrate)	19188	Mucopeptide Biosynthesis
6	F6	Tranexamic Acid	19193	Plasmin-Fibrin Interaction
6	F7	Flibanserin	19203	
6	F8	Naftifine (hydrochloride)	19234	Ergosterol Biosynthesis
6	F9	Fenofibric Acid	19262	PPAR α
6	F10	Primidone	19277	
6	F11	Ceritinib	19374	ALK
6	F12	<i>Unused</i>		
6	G1	<i>Unused</i>		
6	G2	Lenvatinib	19375	VEGFR2, VEGFR3
6	G3	Repaglinide	19387	K _{ir} 6 Channels
6	G4	Clarithromycin	19455	23S rRNA
6	G5	Griseofulvin	19461	Tubulin
6	G6	Capreomycin (sulfate)	19469	
6	G7	Ethacrynic Acid	19536	Na ⁺ /K ⁺ /2Cl ⁻ Symporter
6	G8	Cyproheptadine (hydrochloride hydrate)	19551	5-HT ₂ , Histamine H ₁ , Muscarinic Receptors
6	G9	Cobimetinib	19563	MEK1
6	G10	Podophyllotoxin	19575	Tubulin
6	G11	Argatroban	19639	Thrombin
6	G12	<i>Unused</i>		
6	H1	<i>Unused</i>		
6	H2	Aliskiren (hemifumarate)	19640	Renin
6	H3	Cetirizine (hydrochloride)	19686	Histamine H ₁ Receptors
6	H4	Tolvaptan	19691	Vasopressin V ₂ Receptors
6	H5	AP26113	19778	ALK
6	H6	Aztreonam	19784	Bacterial Cell Wall Synthesis
6	H7	Pomalidomide	19877	Cereblon
6	H8	Tolbutamide	19888	SUR1/K _{IR} 6.2
6	H9	Acalabrutinib	19899	BTK
6	H10	Aripiprazole	19989	
6	H11	Mirtazapine	19994	
6	H12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 17 of 32

Plate	Well	Contents	Item Number	Primary Target*
7	A1	<i>Unused</i>		
7	A2	Eletriptan	20048	5-HT _{1B} , 5-HT _{1D} Receptors
7	A3	Indacaterol	20070	β ₂ -Adrenergic Receptors
7	A4	Diltiazem (hydrochloride)	20079	Calcium Channels
7	A5	Lidocaine	20081	Ion Influx
7	A6	Terazosin (hydrochloride)	20216	α _{1A} -Adrenergic Receptors
7	A7	Idoxuridine	20222	DNA Replication
7	A8	Econazole (nitrate)	20223	CYP51A1
7	A9	Rotigotine	20246	Dopamine D ₁₋₅ Receptors
7	A10	Nelarabine	20248	DNA Replication
7	A11	Labetalol (hydrochloride)	20249	α-Adrenergic, β-Adrenergic Receptors
7	A12	<i>Unused</i>		
7	B1	<i>Unused</i>		
7	B2	Cholic Acid	20250	
7	B3	Nabumetone	20251	COX-1, COX-2
7	B4	N-acetyl-L-Cysteine	20261	
7	B5	Uridine	20300	Thymidylate Synthase
7	B6	Rifapentine	20307	RNA Polymerase
7	B7	Delafloxacin	20314	DNA Gyrase, Topoisomerase IV
7	B8	<i>Unused</i>		
7	B9	Flutamide	20359	Androgen Receptors
7	B10	Betamethasone	20363	Glucocorticoid Receptors
7	B11	Cefadroxil (hydrate)	20365	Penicillin-binding Proteins
7	B12	<i>Unused</i>		
7	C1	<i>Unused</i>		
7	C2	Sulfacetamide (sodium salt)	20377	DHPS
7	C3	Isoniazid	20378	Mycolic Acid Biosynthesis
7	C4	Tetrabenazine	20380	VMAT2
7	C5	Levofloxacin	20382	DNA Gyrase, Topoisomerase IV
7	C6	<i>Unused</i>		
7	C7	Deflazacort	20386	Glucocorticoid Receptors
7	C8	Deferiprone	20387	Iron
7	C9	Flecainide (acetate)	20388	I _{Na} , I _{Kr}
7	C10	Ramelteon	20389	MT ₁ , MT ₂ Receptors
7	C11	Praziquantel	20508	
7	C12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 18 of 32

Plate	Well	Contents	Item Number	Primary Target*
7	D1	<i>Unused</i>		
7	D2	Folic Acid	20515	
7	D3	Ertapenem (sodium salt)	20523	Penicillin-binding Proteins
7	D4	Prednisone	20677	Glucocorticoid Receptors
7	D5	Pentamidine	20679	
7	D6	Acidinium (bromide)	20699	Muscarinic M ₁₋₅ Receptors
7	D7	Fluticasone Propionate	20703	Glucocorticoid Receptors
7	D8	Pyridoxine	20706	
7	D9	Perphenazine	20735	Various CNS Receptors
7	D10	Hydrocortisone	20739	Glucocorticoid, Mineralocorticoid Receptors
7	D11	Risedronate (sodium salt)	20740	Osteoclasts
7	D12	<i>Unused</i>		
7	E1	<i>Unused</i>		
7	E2	Anidulafungin	20753	Glucan Synthase
7	E3	Deoxycholic Acid	20756	
7	E4	Loxapine (succinate)	20760	Various CNS Receptors
7	E5	Piperacillin (sodium salt)	20766	Penicillin-binding Proteins
7	E6	Tolcapone	20768	Catechol O-methyltransferase
7	E7	Etidronate (sodium salt)	20769	Hydroxyapatite, Osteoclasts
7	E8	Perindopril	20796	ACE
7	E9	Etodolac	20833	COX-1, COX-2
7	E10	Acitretin	20853	CRABP-I, CRABP-II
7	E11	Toremifene (citrate)	20854	ER α , ER β
7	E12	<i>Unused</i>		
7	F1	<i>Unused</i>		
7	F2	Prednisolone	20866	Glucocorticoid Receptors
7	F3	Azelastine (hydrochloride)	20873	Histamine H ₁ Receptors
7	F4	Flurandrenolide	20933	EGFR, Glucocorticoid Receptors
7	F5	Norethindrone	20941	Progesterone Receptors
7	F6	Etravirine	20946	Reverse Transcriptase
7	F7	Brompheniramine (maleate)	20996	Histamine H ₁ Receptors
7	F8	Nisoldipine	20998	Calcium Channels
7	F9	Salbutamol	21003	β_2 -Adrenergic Receptors
7	F10	Levobunolol (hydrochloride)	21005	β -Adrenergic Receptors
7	F11	Nafcillin (sodium salt hydrate)	21008	Penicillin-binding Proteins
7	F12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 19 of 32

Plate	Well	Contents	Item Number	Primary Target*
7	G1	<i>Unused</i>		
7	G2	Tipranavir	21028	HIV Protease
7	G3	Gemifloxacin (mesylate)	21047	DNA Gyrase, Topoisomerase IV
7	G4	Vitamin K ₁	21051	Prothrombin, Factor VII, Factor IX, Factor X, Protein C, Protein S, Protein Z
7	G5	Etonogestrel	21062	Progesterone Receptors
7	G6	N-Butyldeoxynojirimycin (hydrochloride)	21065	β -glucosidase II, Ceramide-specific Glucosyltransferase
7	G7	Bacitracin (zinc)	21212	Bacterial Cell Wall Synthesis
7	G8	Acetazolamide	21218	Carbonic Anhydrase
7	G9	Adenosine	21232	Purine Receptors
7	G10	Secnidazole	21236	
7	G11	Mesna	21238	
7	G12	<i>Unused</i>		
7	H1	<i>Unused</i>		
7	H2	Granisetron (hydrochloride)	21239	5-HT ₃ Receptors
7	H3	Triamterene	21242	ENaC
7	H4	(\pm)-Epinephrine (hydrochloride)	21245	α_{1A} -Adrenergic, β_1 -Adrenergic, β_2 -Adrenergic Receptors
7	H5	Guaifenesin	21250	MUC5AC
7	H6	Clobetasol Propionate	21251	Glucocorticoid Receptors
7	H7	Chlorpheniramine (maleate)	21253	Histamine H ₁ Receptors
7	H8	Moexipril (hydrochloride)	21255	ACE
7	H9	Dienogest	21257	Glucocorticoid, Mineralocorticoid Receptors
7	H10	Vinorelbine (tartrate)	21262	Tubulin
7	H11	Enasidenib	21277	IDH2 R140, IDH2 R172
7	H12	<i>Unused</i>		
8	A1	<i>Unused</i>		
8	A2	Tropium (chloride)	21279	Muscarinic M ₁₋₅ Receptors
8	A3	Velpatasvir	21280	NS5A
8	A4	Alcaftadine	21290	Histamine H ₁ Receptors
8	A5	Hydrochlorothiazide	21304	SLC12A3
8	A6	Indapamide	21308	K _v 1.1 Channels
8	A7	Bendroflumethiazide	21311	Na ⁺ /Cl ⁺ Cotransporter
8	A8	Torasemide	21312	Na ⁺ /Cl ⁺ Cotransporter
8	A9	Pantoprazole	21345	H ⁺ /K ⁺ ATPase
8	A10	Ibandronate (sodium hydrate)	21361	Osteoclasts
8	A11	Amantadine (hydrochloride)	21364	NMDA Receptors
8	A12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 20 of 32

Plate	Well	Contents	Item Number	Primary Target*
8	B1	<i>Unused</i>		
8	B2	Mometasone Furoate	21365	Glucocorticoid Receptors
8	B3	Trifluorothymidine	21366	Thymidylate Synthase
8	B4	Desonide	21368	Glucocorticoid Receptors
8	B5	Anagrelide (hydrochloride)	21411	PDE3
8	B6	(-)-Ropivacaine (hydrochloride hydrate)	21422	Sodium Channels
8	B7	Quinapril (hydrochloride)	21439	ACE
8	B8	AN2728	21455	PDE4
8	B9	AHU377 (calcium salt)	21473	Neprilysin
8	B10	VX-661	21480	CFTR
8	B11	Eliglustat (hemitartrate)	21487	Glucosylceramide Synthase
8	B12	<i>Unused</i>		
8	C1	<i>Unused</i>		
8	C2	L-Carnitine	21489	
8	C3	Ertugliflozin	21507	SGLT2
8	C4	Cefdinir	21521	β -Lactamase
8	C5	Retapamulin	21528	Bacterial 50S Ribosomal Subunit
8	C6	Vilazodone (hydrochloride)	21547	5-HT _{1A} Receptors, SERT
8	C7	Rilpivirine	21559	Reverse Transcriptase
8	C8	LY2835219	21560	Cdk4, Cdk6
8	C9	Betrixaban	21563	Factor Xa
8	C10	Afatinib (maleate)	21567	EGFR, HER2
8	C11	Eptifibatide (acetate)	21578	GPIIb, GPIIIa
8	C12	<i>Unused</i>		
8	D1	<i>Unused</i>		
8	D2	Fosamprenavir (calcium salt)	21609	HIV-1 Protease
8	D3	Mycophenolic Acid	21716	IMPDH
8	D4	Megestrol Acetate	21749	Progesterone Receptors
8	D5	Tinidazole	21755	DNA (Oxidative Strand Breaks)
8	D6	Demeclocycline (hydrochloride)	21759	Bacterial 30S Ribosomal Subunit
8	D7	Olodaterol	21794	β_2 -Adrenergic Receptors
8	D8	L(-)- α -Methyldopa	21814	DOPA Decarboxylase
8	D9	Doxylamine (succinate)	21826	Histamine H ₁ Receptors
8	D10	Propafenone (hydrochloride)	21871	Myocardial Membranes
8	D11	Fosinopril (sodium salt)	21899	ACE
8	D12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 21 of 32

Plate	Well	Contents	Item Number	Primary Target*
8	E1	<i>Unused</i>		
8	E2	Dopamine (hydrochloride)	21992	Dopamine Receptors
8	E3	Etomidate	21993	GABA Receptors
8	E4	Buspirone (hydrochloride)	22028	α -Adrenergic Receptors
8	E5	Cangrelor (sodium salt)	22086	P2Y ₁₂ Receptors
8	E6	Simeprevir (sodium salt)	22144	HCV NS3/4A
8	E7	Dolutegravir	22191	HIV Integrase
8	E8	Cysteamine (hydrochloride)	22193	
8	E9	D-Cycloserine	22194	NMDA Receptors
8	E10	Sertaconazole (nitrate)	22232	Sterol 14 α Demethylase
8	E11	Dolasetron	22234	5-HT ₃ Receptors
8	E12	<i>Unused</i>		
8	F1	<i>Unused</i>		
8	F2	Aminophylline	22235	PDE
8	F3	Atomoxetine (hydrochloride)	22248	NET
8	F4	Cabazitaxel	22262	Tubulin
8	F5	Pioglitazone (hydrochloride)	22263	PPAR γ
8	F6	Leuprorelin (acetate)	22275	GnRH Receptor
8	F7	Escitalopram	22405	SERT
8	F8	Alosetron (hydrochloride)	22434	5-HT ₃ Receptors
8	F9	Esmolol (hydrochloride)	22581	β ₁ -Adrenergic Receptors
8	F10	Lifitegrast	22588	LFA-1-ICAM-1 Interaction
8	F11	Avibactam (sodium salt)	22825	β -Lactamase
8	F12	<i>Unused</i>		
8	G1	<i>Unused</i>		
8	G2	Ofloxacin	22891	DNA Gyrase
8	G3	Brexiprazole	22906	5-HT _{1A} , Dopamine D ₂ , D ₃ Receptors
8	G4	Penciclovir	22918	DNA Polymerase
8	G5	Tiagabine (hydrochloride)	22926	GAT-1
8	G6	Malathion	22998	AChE
8	G7	Acetyl- β -methylcholine (chloride)	23092	Muscarinic Receptors
8	G8	AN-2690	23101	tRNA Aminoacylation
8	G9	Ruxolitinib (phosphate)	23215	JAK1, JAK2
8	G10	Disopyramide	23227	Sodium Channels
8	G11	CCK Octapeptide (sulfated)	23371	
8	G12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 22 of 32

Plate	Well	Contents	Item Number	Primary Target*
8	H1	<i>Unused</i>		
8	H2	D-Mannitol	23372	
8	H3	Neostigmine (methyl sulfate)	23501	AChE
8	H4	Nitrofurantoin	23510	Bacterial Ribosomal Proteins
8	H5	Cilastatin	23511	Dipeptidase
8	H6	N-Carbamyl-L-Glutamic Acid	23512	CPS1
8	H7	Tolazamide	23545	SUR1/K _{IR} 6.2
8	H8	Tasimelteon	23546	MT ₁ , MT ₂ Receptors
8	H9	Butoconazole (nitrate)	23554	Sterol 14 α Demethylase
8	H10	Fluphenazine (hydrochloride)	23555	Dopamine D ₁ , D ₂ Receptors
8	H11	Tazarotene	23559	
8	H12	<i>Unused</i>		
9	A1	<i>Unused</i>		
9	A2	Proprantheline (bromide)	23609	Muscarinic Receptors, nAChRs
9	A3	Mivacurium (chloride)	23610	nAChRs
9	A4	Metaproterenol (hemisulfate)	23611	β_2 -Adrenergic Receptors
9	A5	Hydroflumethiazide	23612	Carbonic Anhydrase, Electrolyte Reabsorption
9	A6	Sulfamethoxazole	23613	DHPS
9	A7	Cefotetan (sodium salt)	23624	Bacterial Cell Wall Synthesis
9	A8	Isocarboxazid	23625	MAO
9	A9	Cefaclor	23626	
9	A10	Cefuroxime (sodium salt)	23627	Bacterial Cell Wall Synthesis
9	A11	Fosfomycin Trometamol	23632	MurA
9	A12	<i>Unused</i>		
9	B1	<i>Unused</i>		
9	B2	Cefepime (hydrochloride hydrate)	23633	Bacterial Cell Wall Synthesis
9	B3	<i>Unused</i>		
9	B4	Penicillin V (potassium salt)	23635	Bacterial Cell Wall Synthesis
9	B5	Thiothixene	23649	Dopamine D ₂ Receptors
9	B6	Mefenamic Acid	23650	COX-2
9	B7	Desogestrel	23651	Progesterone Receptors
9	B8	Eslicarbazepine Acetate	23656	Sodium Channels
9	B9	Ulipristal Acetate	23657	Progesterone Receptors
9	B10	Dichlorphenamide	23658	Carbonic Anhydrase
9	B11	Nebivolol (hydrochloride)	23660	β_1 -Adrenergic Receptors
9	B12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Plate	Well	Contents	Item Number	Primary Target*
9	C1	<i>Unused</i>		
9	C2	Olsalazine (sodium salt)	23661	COX, LTB ₄ , Xanthine Oxidase
9	C3	Altretamine	23662	GPX4
9	C4	Histrelin (acetate)	23663	GnRH Receptor
9	C5	Medroxyprogesterone 17-acetate	23664	Progesterone Receptors
9	C6	Mefloquine (hydrochloride)	23665	
9	C7	Pralidoxime (chloride)	23666	AChE
9	C8	Prednicarbate	23668	
9	C9	Ambrisentan	23669	ET _A Receptors
9	C10	6-Mercaptopurine (hydrate)	23675	DNA Synthesis
9	C11	Lesinurad	23688	URAT1
9	C12	<i>Unused</i>		
9	D1	<i>Unused</i>		
9	D2	Pralatrexate	23689	DHFR
9	D3	Besifloxacin (hydrochloride)	23692	Bacterial DNA Gyrase, Topoisomerase IV
9	D4	Bendamustine (hydrochloride)	23693	DNA (Alkylating Agent)
9	D5	Vortioxetine (hydrobromide)	23694	5-HT _{1A} , 5-HT _{1B} , 5-HT _{3A} , 5-HT ₇ Receptors, SERT
9	D6	TAK-599	23696	PBP2
9	D7	Saxagliptin (hydrochloride)	23697	DPP-4
9	D8	Rocuronium (bromide)	23698	nAChRs
9	D9	<i>Unused</i>		
9	D10	Nepafenac	23700	COX-1, COX-2
9	D11	L-Arginine (hydrochloride)	23703	Nitric Oxide
9	D12	<i>Unused</i>		
9	E1	<i>Unused</i>		
9	E2	Albendazole	23705	Tubulin
9	E3	Ethambutol (hydrochloride)	23713	
9	E4	Ethionamide	23714	InhA
9	E5	Didanosine	23715	Viral DNA Reverse Transcriptase
9	E6	L-Glutamine	23716	
9	E7	Mepenzolate (bromide)	23717	Muscarinic M ₂₋₃ Receptors
9	E8	Methimazole	23718	Thyroid Hormone Synthesis
9	E9	Sulfadiazine	23719	DHPS
9	E10	Tiopronin	23720	Cystine
9	E11	Bepotastine (besylate)	23721	Histamine H ₁ Receptors
9	E12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 24 of 32

Plate	Well	Contents	Item Number	Primary Target*
9	F1	<i>Unused</i>		
9	F2	Sulfanilamide	23723	DHPS
9	F3	6-Aminocaproic Acid	23724	Fibrinokinase, Streptokinase, Urokinase
9	F4	Hydroxyurea	23725	
9	F5	4-Aminohippuric Acid	23726	
9	F6	Conivaptan (hydrochloride)	23728	Vasopressin V _{1A} , V ₂ Receptors
9	F7	Tedizolid Phosphate	23729	Bacterial 50S Ribosomal Subunit
9	F8	Daclatasvir	23730	NS5A
9	F9	Naloxegol (oxalate)	23731	μ-Opioid Receptors
9	F10	Ixabepilone	23732	Tubulin
9	F11	D-(+)-Glucose	23733	Energy Metabolism
9	F12	<i>Unused</i>		
9	G1	<i>Unused</i>		
9	G2	Dapsone	23743	
9	G3	Chlorambucil	23744	DNA (Alkylating Agent)
9	G4	<i>Unused</i>		
9	G5	Iohexol	23753	
9	G6	Naratriptan (hydrochloride)	23754	5-HT _{1B} , 5-HT _{1D} Receptors
9	G7	Ospemifene	23755	Estrogen Receptors
9	G8	Octreotide (acetate)	23757	SST ₂ , SST ₃ , SST ₅ Receptors
9	G9	Carteolol (hydrochloride)	23758	β-Adrenergic Receptors
9	G10	Theophylline	23760	Adenosine A ₁ , A ₂ Receptors, PDE
9	G11	Bromfenac (sodium salt)	23763	COX-2
9	G12	<i>Unused</i>		
9	H1	<i>Unused</i>		
9	H2	Peramivir	23765	Neuraminidase
9	H3	Alogliptin (benzoate salt)	23768	DPP-4
9	H4	Dicloxacillin (sodium salt hydrate)	23770	Bacterial Cell Wall Synthesis
9	H5	Frovatriptan (succinate hydrate)	23771	5-HT _{1B} , 5-HT _{1D} Receptors
9	H6	Naphazoline (hydrochloride)	23774	α-Adrenergic Receptors
9	H7	Biperiden (hydrochloride)	23775	Muscarinic Receptors
9	H8	Difluprednate	23776	Glucocorticoid Receptors
9	H9	Fesoterodine (fumarate)	23777	Muscarinic Receptors
9	H10	Pancuronium (bromide hydrate)	23778	nAChRs (Muscle-type)
9	H11	Droxidopa	23779	
9	H12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 25 of 32

Plate	Well	Contents	Item Number	Primary Target*
10	A1	<i>Unused</i>		
10	A2	Halcinonide	23780	
10	A3	Halobetasol Propionate	23781	
10	A4	Mexiletine (hydrochloride)	23782	Sodium Channels
10	A5	Carbidopa (hydrate)	23783	Dopamine Decarboxylase
10	A6	Docosanol	23793	Viral-Host Cell Fusion
10	A7	Doxapram (hydrochloride hydrate)	23794	Carotid Chemoreceptors
10	A8	Lactulose	23795	
10	A9	Retinyl Palmitate	23796	
10	A10	Butenafine (hydrochloride)	23797	Ergosterol Biosynthesis
10	A11	Cortisone Acetate	23798	Glucocorticoid Receptors
10	A12	<i>Unused</i>		
10	B1	<i>Unused</i>		
10	B2	Methyclothiazide	23799	
10	B3	Sulconazole (nitrate)	23800	CYP51A1
10	B4	L-Valacyclovir (hydrochloride)	23801	Viral DNA Polymerase
10	B5	Atovaquone	23802	Complex III
10	B6	Fluocinonide	23803	Glucocorticoid Receptors
10	B7	Benazepril (hydrochloride)	23804	ACE
10	B8	Azilsartan Medoxomil	23805	AT ₂ Receptors
10	B9	Estropipate	23806	OATP1B1
10	B10	Diflorasone Diacetate	23808	Glucocorticoid Receptors
10	B11	Netupitant	23809	NK1
10	B12	<i>Unused</i>		
10	C1	<i>Unused</i>		
10	C2	Mestranol	23822	Estrogen Receptors
10	C3	Oxybutynin (hydrochloride)	23825	Muscarinic Receptors
10	C4	Oxymetazoline (hydrochloride)	23826	5-HT ₁ , α ₁ -Adrenergic, α ₂ -Adrenergic Receptors
10	C5	Bisoprolol (hemifumarate)	23827	β ₁ -Adrenergic Receptors
10	C6	Famotidine	23828	Histamine H ₂ Receptors
10	C7	Acetylcholine (chloride)	23829	Muscarinic Receptors, nAChRs
10	C8	Bethanechol (chloride)	23830	Muscarinic Receptors
10	C9	Pyridostigmine (bromide)	23831	Cholinesterase
10	C10	Methylhomatropine (bromide)	23832	Muscarinic Receptors
10	C11	<i>Unused</i>		
10	C12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 26 of 32

Plate	Well	Contents	Item Number	Primary Target*
10	D1	<i>Unused</i>		
10	D2	Famciclovir	23834	DNA Polymerase
10	D3	Fenopropfen (calcium salt hydrate)	23835	NSAID
10	D4	Avobenzone	23836	UVA
10	D5	Milnacipran (hydrochloride)	23837	NET, SERT
10	D6	Rimantadine (hydrochloride)	23838	
10	D7	Efinaconazole	23839	Ergosterol Biosynthesis
10	D8	Cefprozil	23840	Bacterial Cell Wall Synthesis
10	D9	Molindone (hydrochloride)	23841	Dopamine Receptors
10	D10	Flunisolide	23842	Glucocorticoid Receptors
10	D11	Dihydroergotamine (mesylate)	23847	5-HT _{1B} , 5-HT _{1D} Receptors
10	D12	<i>Unused</i>		
10	E1	<i>Unused</i>		
10	E2	Methylscopolamine (bromide)	23862	Muscarinic Receptors
10	E3	Dicyclomine (hydrochloride)	23867	Muscarinic M ₁₋₂ Receptors
10	E4	Ropinirole (hydrochloride)	23871	Dopamine D ₂ Receptors
10	E5	Fosphenytoin (sodium salt)	23889	Na _v Channels
10	E6	Trimethadione	23893	
10	E7	Amrinone	23898	PDE3
10	E8	Acamprosate (calcium salt)	23899	NMDA Receptors
10	E9	Acetohexamide	23900	SUR1/K _{IR} 6.2
10	E10	Alclometasone Dipropionate	23901	Glucocorticoid Receptors
10	E11	Alvimopan	23902	μ-Opioid Receptors
10	E12	<i>Unused</i>		
10	F1	<i>Unused</i>		
10	F2	Amcinonide	23903	Glucocorticoid Receptors
10	F3	Apraclonidine (hydrochloride)	23904	α ₂ -Adrenergic Receptors
10	F4	Cetrorelix (acetate)	23910	GnRH Receptor
10	F5	Cabergoline	23934	Dopamine Receptors
10	F6	Arformoterol (tartrate)	23935	β ₂ -Adrenergic Receptors
10	F7	Benzonatate	23936	Na _v Channels
10	F8	Carbinoxamine (maleate)	23937	Histamine H ₁ Receptors
10	F9	Metaraminol (tartrate)	23938	β ₁ -Adrenergic Receptors
10	F10	Crotamiton	23943	TRPV4
10	F11	Cyclopentolate (hydrochloride)	23944	Muscarinic M ₁₋₃ Receptors
10	F12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 27 of 32

Plate	Well	Contents	Item Number	Primary Target*
10	G1	<i>Unused</i>		
10	G2	Desoximetasone	23945	Glucocorticoid Receptors
10	G3	Emedastine (fumarate)	23946	Histamine H ₁ Receptors
10	G4	Ethosuximide	23947	
10	G5	Hexachlorophene	23948	D-Lactate Dehydrogenase
10	G6	Hexaminolevulinate (hydrochloride)	23949	
10	G7	Isavuconazonium (sulfate)	23950	Ergosterol Biosynthesis
10	G8	Levomilnacipran (hydrochloride)	23951	NET, SERT
10	G9	γ-Lindane	23952	GABAA Receptor/Chloride Ionophore Complex
10	G10	Nilutamide	23953	Androgen Receptors
10	G11	Oxacillin (sodium salt)	23954	Bacterial Cell Wall Synthesis
10	G12	<i>Unused</i>		
10	H1	<i>Unused</i>		
10	H2	D-Penicillamine	23955	Copper
10	H3	Phenelzine (sulfate)	23956	MAO
10	H4	Quinine	23958	<i>P. falciparum</i> Nucleic Acid Synthesis, Protein Synthesis, Glycolysis
10	H5	Terconazole	23959	Ergosterol Biosynthesis
10	H6	Azelaic Acid	23977	Multiple Enzymes
10	H7	Hydroxyprogesterone Caproate	23978	Glucocorticoid Receptors
10	H8	Acetohydroxamic Acid	23979	Urease
10	H9	Isoflurane	23989	GABA _A Receptors
10	H10	Isosorbide Dinitrate	23990	Nitric Oxide
10	H11	Salbutamol (hydrochloride)	23991	β ₂ -Adrenergic Receptors
10	H12	<i>Unused</i>		
11	A1	<i>Unused</i>		
11	A2	(R)-Cetirizine (hydrochloride)	23992	Histamine H ₁ Receptors
11	A3	Levonordefrin	23993	α ₂ -Adrenergic Receptors
11	A4	Lodoxamide	23994	GPR35, Mast Cells
11	A5	Mafenide	23995	
11	A6	Sevoflurane	23996	GABA _A , Glycine Receptors
11	A7	Trientine (hydrochloride)	24000	Copper
11	A8	Methyldopate	24002	Dopamine Decarboxylase
11	A9	<i>Unused</i>		
11	A10	Metipranolol (hydrochloride)	24003	β ₂ -Adrenergic Receptors
11	A11	Methenamine Hippurate	24004	
11	A12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 28 of 32

Plate	Well	Contents	Item Number	Primary Target*
11	B1	<i>Unused</i>		
11	B2	Nedocromil (sodium salt)	24005	Mast Cells
11	B3	Oxtriphylline	24006	Adenosine A ₁ , A ₂ Receptors, PDE
11	B4	Tamsulosin (hydrochloride)	24020	α ₁ -Adrenergic Receptors
11	B5	Acrivastine	24021	Histamine H ₁ Receptors
11	B6	Ethotoin	24022	Sodium Ion Influx
11	B7	Cariprazine (hydrochloride)	24025	Various CNS Receptors
11	B8	Delavirdine (mesylate)	24026	HIV-1 Reverse Transcriptase
11	B9	Dexchlorpheniramine (maleate)	24027	Histamine H ₁ Receptors
11	B10	Dipivefrin (hydrochloride)	24028	α _{1A} -Adrenergic, β ₁ -Adrenergic, β ₂ -Adrenergic Receptors
11	B11	Echothiophate (iodide)	24029	AChE
11	B12	<i>Unused</i>		
11	C1	<i>Unused</i>		
11	C2	Elbasvir	24031	NS5A
11	C3	Fludrocortisone Acetate	24032	Mineralocorticoid Receptors
11	C4	<i>Unused</i>		
11	C5	Fosaprepitant dimeglumine	24034	NK1
11	C6	Phenytoin	24037	Na _v Channels
11	C7	Degarelix (acetate)	24069	GnRH Receptor
11	C8	Nafarelin (acetate)	24070	GnRH Receptor
11	C9	Methylnaltrexone (bromide)	24072	μ-Opioid Receptors
11	C10	Linaclotide	24085	Guanylate Cyclase C Receptor
11	C11	Oritavancin (phosphate)	24091	
11	C12	<i>Unused</i>		
11	D1	<i>Unused</i>		
11	D2	Ergoloid (mesylates)	24095	
11	D3	Sugammadex (sodium salt)	24096	
11	D4	Ganirelix (acetate)	24098	GnRH Receptor
11	D5	Hydroxocobalamin	24099	
11	D6	Trimethobenzamide (hydrochloride)	24145	Dopamine D ₂ Receptors
11	D7	P005672 (hydrochloride)	24685	Bacterial 70S Ribosomal Subunit
11	D8	Miltefosine	63280	CTP:phosphocholine Cytidylyltransferase
11	D9	(±)-Flurbiprofen	70250	COX-1, COX-2
11	D10	Aspirin	70260	COX-1, COX-2
11	D11	5-Aminosalicylic Acid	70265	
11	D12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 29 of 32

Plate	Well	Contents	Item Number	Primary Target*
11	E1	<i>Unused</i>		
11	E2	Indomethacin	70270	COX-1, COX-2
11	E3	(±)-Ibuprofen	70280	COX-1, COX-2
11	E4	(+)-Naproxen	70290	COX-1, COX-2
11	E5	Meclofenamate (sodium salt)	70550	COX-1, COX-2
11	E6	Diclofenac (sodium salt)	70680	COX-1, COX-2
11	E7	Ketorolac (tromethamine salt)	70690	COX-1, COX-2
11	E8	Rosiglitazone	71740	PPAR γ
11	E9	Calcitriol	71820	Vitamin D Receptor
11	E10	(6R)-5,6,7,8-tetrahydro-L-Biopterin (hydrochloride)	81880	Phenylalanine Hydroxylase, Tryptophan Hydroxylase, Tyrosine Hydroxylase
11	E11	Oleoyl Ethanolamide	90265	PPAR α
11	E12	<i>Unused</i>		
11	F1	<i>Unused</i>		
11	F2	Capsaicin	92350	TRPV1
11	F3	Fluorometholone	9000157	Glucocorticoid Receptors
11	F4	Spirolactone	9000324	Aldosterone, Androgen, Mineralocorticoid Receptors
11	F5	(±)-Asenapine	9000496	Various CNS Receptors
11	F6	Dronedarone (hydrochloride)	9000543	I $_K$
11	F7	Riociguat	9000554	sGC
11	F8	Lurasidone	9000570	Various CNS Receptors
11	F9	<i>Unused</i>		
11	F10	25-hydroxy Vitamin D ₃	9000683	
11	F11	Colchicine	9000760	Tubulin
11	F12	<i>Unused</i>		
11	G1	<i>Unused</i>		
11	G2	Vigabatrin	9000976	GABA-T
11	G3	Levetiracetam	9001820	SV2A
11	G4	4-hydroxy Amphetamine (hydrochloride)	9001830	TAAR1
11	G5	PF-299804	9001879	EGFR, HER2, HER4
11	G6	(±)-O-Desmethyl Venlafaxine	9001881	NET, SERT
11	G7	Saquinavir (mesylate)	9001893	HIV Protease
11	G8	5-Aminolevulinic Acid (hydrochloride)	9001902	
11	G9	Cephalexin	9002009	PBP3
11	G10	Metronidazole	9002409	
11	G11	Sulindac	10004386	COX-1, COX-2
11	G12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 30 of 32

Plate	Well	Contents	Item Number	Primary Target*
11	H1	<i>Unused</i>		
11	H2	Fenofibrate	10005368	PPAR α
11	H3	Orlistat	10005426	Lipases
11	H4	Taflopust	10005440	FP Receptors
11	H5	17 β -Estradiol	10006315	ER α , ER β
11	H6	Mifepristone	10006317	Androgen, Glucocorticoid, Progesterone Receptors
11	H7	Levonorgestrel	10006318	Androgen, Progesterone Receptors
11	H8	Ethynyl Estradiol	10006486	ER α , ER β
11	H9	Losartan (potassium salt)	10006594	AT $_1$ Receptors
11	H10	(\pm)-Ketoprofen	10006661	COX-1, COX-2
11	H11	Zileuton	10006967	5-LO
11	H12	<i>Unused</i>		
12	A1	<i>Unused</i>		
12	A2	FK-506	10007965	FKBP12
12	A3	Zafirlukast	10008282	CysLT $_1$
12	A4	Montelukast (sodium salt)	10008318	CysLT $_1$
12	A5	Gabapentin	10008346	
12	A6	Celecoxib	10008672	COX-2
12	A7	Bortezomib	10008822	20S Proteasome
12	A8	Eicosapentaenoic Acid ethyl ester	10008884	
12	A9	Calcipotriol (hydrate)	10009599	VDR
12	A10	Lomitapide	10009610	MTTP
12	A11	Sorafenib	10009644	Multiple Kinases
12	A12	<i>Unused</i>		
12	B1	<i>Unused</i>		
12	B2	SAHA	10009929	HDAC
12	B3	Fluvastatin (sodium salt hydrate)	10010337	HMG-CoA Reductase
12	B4	Lovastatin	10010338	HMG-CoA Reductase
12	B5	Pravastatin (sodium salt)	10010343	HMG-CoA Reductase
12	B6	Simvastatin	10010344	HMG-CoA Reductase
12	B7	NS 304	10010411	IP Receptors
12	B8	Nilotinib	10010422	Bcr-Abl
12	B9	Lubiprostone	10010487	CIC-2
12	B10	Tranylcypromine (hydrochloride)	10010494	LSD1, MAO-A, MAO-B
12	B11	Fulvestrant	10011269	ER α
12	B12	<i>Unused</i>		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 31 of 32

Plate	Well	Contents	Item Number	Primary Target*
12	C1	Unused		
12	C2	Plerixafor (hydrochloride hydrate)	10011332	CXCR4
12	C3	Terbinafine (hydrochloride)	10011619	Squalene Epoxidase
12	C4	Raloxifene (hydrochloride)	10011620	ER α , Er β , K ν 4.3
12	C5	Allopurinol	10012597	Xanthine Oxidoreductase
12	C6	Nefazodone (hydrochloride)	10012642	5-HT $_{2A}$, α_2 -Adrenergic, Dopamine D $_2$ Receptors
12	C7	Unused		
12	C8	Unused		
12	C9	Unused		
12	C10	Unused		
12	C11	Unused		
12	C12	Unused		
12	D1	Unused		
12	D2	Unused		
12	D3	Unused		
12	D4	Unused		
12	D5	Unused		
12	D6	Unused		
12	D7	Unused		
12	D8	Unused		
12	D9	Unused		
12	D10	Unused		
12	D11	Unused		
12	D12	Unused		
12	E1	Unused		
12	E2	Unused		
12	E3	Unused		
12	E4	Unused		
12	E5	Unused		
12	E6	Unused		
12	E7	Unused		
12	E8	Unused		
12	E9	Unused		
12	E10	Unused		
12	E11	Unused		
12	E12	Unused		

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 32 of 32

Plate	Well	Contents	Item Number	Primary Target*
12	F1	Unused		
12	F2	Unused		
12	F3	Unused		
12	F4	Unused		
12	F5	Unused		
12	F6	Unused		
12	F7	Unused		
12	F8	Unused		
12	F9	Unused		
12	F10	Unused		
12	F11	Unused		
12	F12	Unused		
12	G1	Unused		
12	G2	Unused		
12	G3	Unused		
12	G4	Unused		
12	G5	Unused		
12	G6	Unused		
12	G7	Unused		
12	G8	Unused		
12	G9	Unused		
12	G10	Unused		
12	G11	Unused		
12	G12	Unused		
12	H1	Unused		
12	H2	Unused		
12	H3	Unused		
12	H4	Unused		
12	H5	Unused		
12	H6	Unused		
12	H7	Unused		
12	H8	Unused		
12	H9	Unused		
12	H10	Unused		
12	H11	Unused		
12	H12	Unused		

*Primary Target is based on available IC_{50} and/or K_i values cited in the literature

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM