

PRODUCT INFORMATION



Cellular Metabolism Screening Library

Item No. 33705 • Batch No. 0609421

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

Page 1 of 6

Plate	Well	Contents	Item Number	Primary Target*
1	A1	<i>Unused</i>		
1	A2	TMS	10038	CYP1A1
1	A3	BEC	10170	Arginase I, II
1	A4	Paclitaxel	10461	DNA Synthesis
1	A5	PK11195	10525	TPSO
1	A6	Olaparib	10621	PARP1, PARP2
1	A7	SC-26196	10792	Δ^6 Desaturase
1	A8	Torin 1	10997	mTOR
1	A9	Piperlongumine	11006	GSTP1
1	A10	17-DMAG	11036	Hsp90
1	A11	STF-31	11173	Glut1
1	A12	<i>Unused</i>		
1	B1	<i>Unused</i>		
1	B2	Actinomycin D	11421	DNA Synthesis
1	B3	ABT-263	11500	Bcl-2, Bcl-xL, Bcl-W
1	B4	GSK1059615	11569	PI3K α
1	B5	Phlorizin	11576	SGLT1, SGLT2
1	B6	Betulinic Acid	11686	TGR5
1	B7	Gemcitabine	11690	Cytidine
1	B8	6-Gingerol	11707	TRPV1, TRPA1
1	B9	Artemisinin	11816	Antimalarial
1	B10	Artesunate	11817	GST
1	B11	6-Shogaol	11901	Mitochondrial Uncoupler
1	B12	<i>Unused</i>		
1	C1	<i>Unused</i>		
1	C2	Dorsomorphin	11967	AMPK
1	C3	(+)-Etomoxir (sodium salt)	11969	CPT1
1	C4	Anastrozole	11987	CYP19A1
1	C5	Rottlerin	12006	Mitochondrial Uncoupler
1	C6	Rosuvastatin	12029	HMGCR
1	C7	CHIR99021	13122	GSK3
1	C8	JZL 184	13158	MAGL
1	C9	FK-866	13287	NAMPT
1	C10	Fatostatin (hydrobromide)	13562	SREBP
1	C11	YM-201636	13576	PIKfyve
1	C12	<i>Unused</i>		

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, 03/01/2024

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



Plate	Well	Contents	Item Number	Primary Target*
1	D1	<i>Unused</i>		
1	D2	OU749	13804	GGT
1	D3	ML-265	13942	PKM2
1	D4	Methotrexate (hydrate)	13960	DHFR
1	D5	Rotenone	13995	Complex I
1	D6	AK-7	14004	SIRT2
1	D7	G6PDi-1	31484	G6PDH
1	D8	Febuxostat	14127	Xanthine Oxidase
1	D9	Fludarabine	14128	Purine Analog
1	D10	Plumbagin	14314	NOX
1	D11	AESBF (hydrochloride)	14321	NOX, Serine Proteases
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	5-Fluorouracil	14416	Thymidylate Synthase
1	E3	Gossypol	14482	LDH, AKR1B1
1	E4	Dorzolamide (hydrochloride)	14616	CAII, CAIV
1	E5	AGI-5198	14624	IDH1 R132H
1	E6	Bumetanide	14630	NKCC1, NKCC2
1	E7	AGI-6780	14639	IDH2 R140Q
1	E8	Lonidamine	14640	HK
1	E9	3-Nitropropionic Acid	14684	Complex II
1	E10	Dimethyl fumarate	14714	GSH
1	E11	TBCA	14831	CK2
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	Pentostatin	14878	Adenosine Deaminase
1	F3	Finasteride	14938	5 α -reductase II
1	F4	Metyrapone	14994	Cortisol
1	F5	Linezolid	15012	Protein Synthesis
1	F6	Miglitol	15014	α -Glucosidase
1	F7	PF-04981517	15019	CYP3A4
1	F8	Torcetrapib	15028	CETP
1	F9	PF-429242 (hydrochloride)	15140	Site-1 Protease
1	F10	HQNO	15159	Complex III
1	F11	FCCP	15218	Mitochondrial Uncoupler
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 6

Plate	Well	Contents	Item Number	Primary Target*
1	G1	<i>Unused</i>		
1	G2	Atglistatin	15284	ATGL
1	G3	Auranofin	15316	TxR1
1	G4	SF1670	15368	PTEN
1	G5	Ranolazine (hydrochloride)	15604	Sodium Channels
1	G6	CGP 37157	15611	Mitochondrial Sodium-Calcium Exchanger
1	G7	6-Thioguanine	15774	Purine Analog
1	G8	WWL113	16181	CES3, CES1F
1	G9	BMS 303141	16239	ACL
1	G10	PR-619	16276	DUBs
1	G11	TG101348	16289	JAK2, THTR2
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	MitoTEMPO (hydrate)	16621	SOD Mimetic
1	H3	U-104	16717	CAIX, CAXII
1	H4	Swainsonine	16860	Lysosomal α -Mannosidase, Golgi α -Mannosidase II
1	H5	DL- α -Difluoromethylornithine (hydrochloride hydrate)	16889	ODC
1	H6	UK 5099	16980	MPC
1	H7	CPI-613	16981	PDH, α -Ketoglutarate Dehydrogenase
1	H8	Toyocamycin (hydrate)	17371	IRE1, PI3K, Rio1 Kinase
1	H9	GSK2656157	17372	PERK
1	H10	Vps34-IN1	17392	VPS34
1	H11	6-diazo-5-oxo-L-nor-Leucine	17580	Glutaminase, CTPS
1	H12	<i>Unused</i>		
2	A1	<i>Unused</i>		
2	A2	Fenretinide	17688	DEGS
2	A3	PFK15	17689	PFKFB3
2	A4	Erastin	17754	System Xc
2	A5	Spautin-1	17769	USP10 and USP13
2	A6	Fomepizole	17851	ADH1
2	A7	Nitisinone	17924	4-HPPD
2	A8	Lometrexol (hydrate)	18049	GARFT
2	A9	Dipyridamole	18189	PDE, ENT
2	A10	CVT-10216	18318	ALDH2
2	A11	GDC-0068	18412	Akt
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



Plate	Well	Contents	Item Number	Primary Target*
2	B1	<i>Unused</i>		
2	B2	S3QEL-2	18556	Complex III
2	B3	Mitotane	18651	CYP11B1
2	B4	2-Acetylphenothiazine	19056	NOX1
2	B5	A-939572	19123	SCD1
2	B6	NK 252	19131	Nrf2
2	B7	VAS2870	19205	NOX
2	B8	AZD 7545	19282	PDK1, PDK3
2	B9	BPTES	19284	GLS
2	B10	CB-5083	19311	p97 ATPase
2	B11	NCT-503	19718	PHGDH
2	B12	<i>Unused</i>		
2	C1	<i>Unused</i>		
2	C2	Adaptaquin	19720	PHD2
2	C3	WZB117	19900	Glut1
2	C4	AZD 3965	19912	MCT1
2	C5	GSK2194069	20022	FASN
2	C6	Copanlisib	20354	PI3K α , β , δ , γ
2	C7	Deferiprone	20387	Iron
2	C8	GSK2837808A	20626	LDHA/B
2	C9	Dicoumarol	20764	NQO1, SAPK
2	C10	Tolcapone	20768	COMT
2	C11	BTB06584	21032	F ₁ F ₀ ATPase
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	NCT-501	21162	ALDH1A1
2	D3	R-59-949	21221	DGK- α
2	D4	Eliglustat (hemitartrate)	21487	GCS
2	D5	1-Deoxynojirimycin	21500	α -Glucosidase
2	D6	(\pm)-NLG-919	21509	IDO Pathway
2	D7	Mycophenolic Acid	21716	IMPDH
2	D8	CJ-13610	21764	5-LO
2	D9	PFK158	22987	PFKFB3
2	D10	ML-210	23282	GPX4
2	D11	Tofogliflozin (hydrate)	23509	SGLT2
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



Plate	Well	Contents	Item Number	Primary Target*
2	E1	<i>Unused</i>		
2	E2	Mefloquine (hydrochloride)	23665	Lysosomes
2	E3	6-Mercaptopurine (hydrate)	23675	Purine Synthesis
2	E4	Orteronel	24191	CYP17A1
2	E5	ARN14988	24284	Acid Ceramidase
2	E6	Brequinar	24445	DHODH
2	E7	Raltitrexed	26079	Thymidylate Synthase
2	E8	CCF-642	27054	PDI
2	E9	Digitoxin	27825	HIF-1 α
2	E10	DS44960156	28366	MTHFD2
2	E11	Bisubstrate Inhibitor 78	28567	NNMT
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	iFSP1	29483	FSP1
2	F3	Verinurad	29667	URAT1
2	F4	Alrestatin	29888	AKR1B1
2	F5	Pantothenate Kinase Inhibitor	31002	PanK
2	F6	PRGL493	32748	ACSL4
2	F7	Myriocin	63150	Serine Palmitoyltransferase
2	F8	Miltefosine	63280	CTP:PCT
2	F9	SC-560	70340	COX-1
2	F10	Ebselen	70530	GPX mimic
2	F11	Chlorogenic Acid	70930	Phenol Natural Product
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	Rosiglitazone	71740	PPAR γ
2	G3	Diphenyleiodonium (chloride)	81050	iNOS, eNOS
2	G4	NS 2028	81600	Soluble Guanylyl Cyclase
2	G5	Oleoyl Ethanolamide	90265	PPAR α
2	G6	Vigabatrin	9000976	GABA-T
2	G7	Orlistat	10005426	DAGL α , Lipases
2	G8	EUK 134	10006329	SOD and Catalase Mimetic
2	G9	Ro 48-8071	10006415	OSC
2	G10	GW 0742	10006798	PPAR δ
2	G11	AS-605240	10007707	PI3K $\alpha,\beta,\delta,\gamma$
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 6

Plate	Well	Contents	Item Number	Primary Target*
2	H1	Unused		
2	H2	Celecoxib	10008672	COX-2
2	H3	Sphingosine Kinase Inhibitor 2	10009222	SPHK1
2	H4	6-Aminonicotinamide	10009315	G6PD
2	H5	(±)-EX-527	10009798	SIRT1
2	H6	BMS309403	10010206	FABP4
2	H7	TMP-153	10010516	ACAT
2	H8	β-cyano-L-Alanine	10010947	CSE
2	H9	Eeyarestatin 1	10012609	ERAD
2	H10	A-922500	10012708	DGAT-1
2	H11	ND-630	23961	ACC1, ACC2
2	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