

PRODUCT INFORMATION



Antiviral Screening Library

Item No. 30390 • Batch No. 0586543

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

Page 1 of 18

Plate	Well	Contents	Item Number	Primary Target*
1	A1	<i>Unused</i>		
1	A2	Fumagillin	11332	Antiparasitic
1	A3	Hypericin	11334	Broad-spectrum Antiviral
1	A4	5'-Ethylnyl-2'-deoxycytidine	11581	HSV and Vaccinia Virus
1	A5	Aloe-emodin	11677	Undefined Antiviral Association
1	A6	Andrographolide	11679	Antiviral
1	A7	Psoralen	11751	Undefined Antiviral Association
1	A8	(-)-Epicatchin Gallate	11808	Undefined Antiviral Association
1	A9	Artemisinin	11816	Undefined Antiviral Association
1	A10	Artesunate	11817	Undefined Antiviral Association
1	A11	Corilagin	11832	Undefined Antiviral Association
1	A12	<i>Unused</i>		
1	B1	<i>Unused</i>		
1	B2	Glabridin	11843	Antiviral
1	B3	Glycyrrhizic Acid (ammonium salt)	11847	Antiviral
1	B4	Kaempferol	11852	Undefined Antiviral Association
1	B5	Licochalcone A	11853	Undefined Antiviral Association
1	B6	AMD 3465 (hydrobromide)	11959	HIV
1	B7	Oxyresveratrol	12028	Undefined Antiviral Association
1	B8	Astragaloside IV	12069	Adenovirus Type 3
1	B9	Daidzin	13202	Undefined Antiviral Association
1	B10	YM-201636	13576	HIV
1	B11	Ritonavir	13872	HIV-1 Protease
1	B12	<i>Unused</i>		
1	C1	<i>Unused</i>		
1	C2	Nervonic Acid	13940	HIV-1 Reverse Transcriptase
1	C3	Dynasore	14062	Undefined Antiviral Association
1	C4	Glycitin	14161	Undefined Antiviral Association
1	C5	AEBSF (hydrochloride)	14321	Undefined Antiviral Association
1	C6	S3I-201	14336	STAT3
1	C7	Honokiol	14597	Undefined Antiviral Association
1	C8	DMXAA	14617	Undefined Antiviral Association
1	C9	Salubrinal	14735	Herpes Virus Replication
1	C10	R-848	14806	TLR7, TLR8
1	C11	Mitoxantrone (hydrochloride)	14842	HIV-1 Integrase
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, 04/21/2021

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 18

Plate	Well	Contents	Item Number	Primary Target*
1	D1	<i>Unused</i>		
1	D2	ZCL 278	14849	Viral Entry Inhibitor, Envelope Viruses
1	D3	Arctigenin	14913	Influenza
1	D4	Oxymatrine	14915	HBV
1	D5	Oltipraz	15118	Undefined Antiviral Association
1	D6	Q-VD-OPH	15260	Antiviral, <i>pan</i> -Caspase Inhibitor
1	D7	Ginsenoside Rb ₁	15319	Undefined Antiviral Association
1	D8	Harringtonine	15361	Undefined Antiviral Association
1	D9	Arctiin	15375	Influenza
1	D10	FR180204	15544	Dengue Virus
1	D11	Loratadine	15625	Undefined Antiviral Association
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	Oseltamivir Acid	15779	Neuraminidase
1	E3	Fialuridine	15867	Viral Thymidine Kinase
1	E4	N-Acetylneuraminic Acid	16091	Undefined Antiviral Association
1	E5	BIO-Acetoxime	16329	Herpes Viruses
1	E6	Ku-55933	16336	HIV-1 Replication
1	E7	2,4-Diacetylphloroglucinol	16345	Antiviral
1	E8	Monensin (sodium salt)	16488	Viral Replication
1	E9	Isoprinosine	16492	Undefined Antiviral Association
1	E10	3'-Sialyllactose (sodium salt)	16617	Antiviral
1	E11	Umifenovir (hydrochloride)	16933	Broad-spectrum Viral, Fusion Inhibitor
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	Casticin	16976	Undefined Antiviral Association
1	F3	GSK650394	17001	Undefined Antiviral Association
1	F4	ML-324	17472	JMJD2E
1	F5	Hinokitiol	17538	Undefined Antiviral Association
1	F6	Letermovir	17556	CMV replication
1	F7	Formononetin	17558	Antiviral
1	F8	Clemizole (hydrochloride)	17695	HCV NS4B
1	F9	L-Cycloserine	17780	Undefined Antiviral Association
1	F10	VP-14637	17805	RSV Fusion
1	F11	Tilorone (hydrochloride)	17868	Undefined Antiviral Association
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 18

Plate	Well	Contents	Item Number	Primary Target*
1	G1	<i>Unused</i>		
1	G2	Elacridar	18128	HIV
1	G3	Vidarabine	18149	DNA polymerase
1	G4	Boceprevir	18379	HCV NS3/4A
1	G5	Golgicide A	18430	Antiviral
1	G6	Patchouli Alcohol	18450	Antiviral
1	G7	Quercetin 3-D-galactoside	18648	Undefined Antiviral Association
1	G8	Pentoxifylline	18720	HIV
1	G9	Gardiquimod	18877	TLR7, HIV-1
1	G10	ESI-09	19130	Undefined Antiviral Association
1	G11	Osthole	19195	Undefined Antiviral Association
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	Chlorcyclizine (hydrochloride)	19239	HCV
1	H3	CAY10704	19240	HCV
1	H4	Ambroxol (hydrochloride)	19377	Influenza
1	H5	Monoacetylphloroglucinol (hydrate)	19451	Antiviral Precursor
1	H6	D,L-1'-Acetoxychavicol Acetate	19649	Viral Replication
1	H7	Birinapant	19699	HBV
1	H8	Baicalin	19843	HIV-1 Reverse Transcriptase
1	H9	Schisandrin A	19849	Undefined Antiviral Association
1	H10	Telaprevir	20054	HCV NS3/4A
1	H11	Etidronate (sodium salt)	20769	Undefined Antiviral Association
1	H12	<i>Unused</i>		
2	A1	<i>Unused</i>		
2	A2	(+)-JQ1	11187	IFN- γ Suppressor
2	A3	Saracatinib	11497	Dengue Virus RNA Replication
2	A4	CP 690,550	11598	Undefined Antiviral Association
2	A5	Betulinic Acid	11686	HIV-1 Replication
2	A6	Oleanolic Acid	11726	Antiviral
2	A7	Atazanavir	11733	HIV-1 Protease
2	A8	Neoandrographolide	11742	Antiviral
2	A9	Brefeldin A	11861	Undefined Antiviral Association
2	A10	Valproic Acid (sodium salt)	13033	Undefined Antiviral Association
2	A11	Cidofovir	13113	Herpes Viruses
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 18

Plate	Well	Contents	Item Number	Primary Target*
2	B1	<i>Unused</i>		
2	B2	Imatinib (mesylate)	13139	Undefined Antiviral Association
2	B3	Panobinostat	13280	HIV
2	B4	Itraconazole	13288	Undefined Antiviral Association
2	B5	Nitazoxanide	13692	Undefined Antiviral Association
2	B6	Ganciclovir	13853	CMV
2	B7	Lopinavir	13854	HIV-1 Protease
2	B8	Tenofovir	13874	HIV Reverse Transcriptase, HBV Polymerase
2	B9	Valganciclovir (hydrochloride)	13875	Viral DNA Synthesis
2	B10	Acyclovir	14160	Viral DNA Polymerase
2	B11	Chloroquine (phosphate)	14194	Corona Virus (COVID-19)
2	B12	<i>Unused</i>		
2	C1	<i>Unused</i>		
2	C2	Pemetrexed (sodium salt hydrate)	14269	Undefined Antiviral Association
2	C3	2-deoxy-D-Glucose	14325	Undefined Antiviral Association
2	C4	Efavirenz	14412	Reverse Transcriptase
2	C5	5-Fluorouracil	14416	Undefined Antiviral Association
2	C6	Maraviroc	14641	CCR5
2	C7	Abacavir (sulfate)	14746	HIV-1 Reverse Transcriptase
2	C8	Shikonin	14751	HIV-1 Replication
2	C9	Imiquimod	14956	TLR7
2	C10	Stavudine	14975	HIV Reverse Transcriptase
2	C11	Probenecid	14981	Undefined Antiviral Association
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	Doxorubicin (hydrochloride)	15007	HIV
2	D3	Trifluoperazine (hydrochloride)	15068	Influenza
2	D4	Nevirapine	15117	Reverse Transcriptase
2	D5	Zanamivir	15123	Neuraminidase
2	D6	Roflumilast	15141	Undefined Antiviral Association
2	D7	Nelfinavir (mesylate)	15144	HIV-1 Protease
2	D8	DBeQ	15318	Undefined Antiviral Association
2	D9	Amprenavir	15369	HIV Protease
2	D10	Zerumbone	15400	Epstein-Barr
2	D11	Zidovudine	15492	Reverse Transcriptase
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 18

Plate	Well	Contents	Item Number	Primary Target*
2	E1	<i>Unused</i>		
2	E2	6-Thioguanine	15774	Undefined Antiviral Association
2	E3	Darunavir	15866	HIV-1 Protease
2	E4	Cytarabine	16069	DNA Polymerase
2	E5	Oseltamivir (phosphate)	16070	Influenza
2	E6	Raltegravir (potassium salt)	16071	HIV-1 Integrase
2	E7	RVX-208	16424	HIV
2	E8	AZ 960	16731	JAK
2	E9	Ribavirin	16757	DNA and RNA Viruses
2	E10	Emtricitabine	16879	Influenza
2	E11	Tenofovir Disoproxil (fumarate)	16922	Prodrug, HIV Reverse Transcriptase, HBV Polymerase Inhibitor
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	Limonin	16932	HIV
2	F3	OG-L002	17471	LSD1
2	F4	Spautin-1	17769	Undefined Antiviral Association
2	F5	Elvitegravir	17798	HIV-1 Integrase
2	F6	Lamivudine	18514	HIV-1, HIV-2, HBV Reverse Transcriptase
2	F7	Chrysophanol	19870	HBV
2	F8	Idoxuridine	20222	HSV-1 DNA Replication
2	F9	N-acetyl-L-Cysteine	20261	Undefined Antiviral Association
2	F10	Deferiprone	20387	HIV
2	F11	Risedronate (sodium salt)	20740	Undefined Antiviral Association
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	Asunaprevir	20835	HCV NS3/4A
2	G3	Nomilin	20836	HIV-1 Replication
2	G4	Cephaeline	20916	Zika Virus
2	G5	Etravirine	20946	Reverse Transcriptase
2	G6	N-Butyldeoxyjirimycin (hydrochloride)	21065	HIV-1, HIV-2
2	G7	NCH-51	21234	HIV-1 Replication
2	G8	NBD-556	21274	HIV-1 Viral Entry
2	G9	Velpatasvir	21280	HCV NS5A
2	G10	Telbivudine	21453	HBV
2	G11	Bicyclol	21551	HBV
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 18

Plate	Well	Contents	Item Number	Primary Target*
2	H1	<i>Unused</i>		
2	H2	Rilpivirine	21559	Reverse Transcriptase
2	H3	(-)- β -Pinene	21577	Infectious Bronchitis Virus Replication
2	H4	Fosamprenavir (calcium salt)	21609	Prodrug, HIV-1 Protease Inhibitor
2	H5	Moroxydine (hydrochloride)	21647	Broad-spectrum Antiviral
2	H6	Mycophenolic Acid	21716	Antiviral
2	H7	Megestrol Acetate	21749	HIV
2	H8	Amentoflavone	21779	Influenza A, Influenza B, HSV-1
2	H9	Grazoprevir	21816	HCV NS3/4A
2	H10	Pirlindole	22060	Coxsackievirus B3 Genome Replication, Enterovirus D68
2	H11	BAY 57-1293	22129	HSV Helicase-Primase
2	H12	<i>Unused</i>		
3	A1	<i>Unused</i>		
3	A2	Simeprevir (sodium salt)	22144	HCV NS3/4A
3	A3	Dolutegravir	22191	HIV Integrase
3	A4	Ginsenoside Rb ₂	22265	Undefined Antiviral Association
3	A5	Mangiferin	22360	HIV-1
3	A6	Procyanidin B1	22411	HCV RNA Replication
3	A7	2'-C-Methylcytidine	22887	Broad-spectrum Antiviral
3	A8	Matrine	22901	HBV
3	A9	YYA-021	22915	HIV Viral Entry
3	A10	Penciclovir	22918	HSV-1, HSV-2, VZV
3	A11	(1R)-(+)-Camphor	23175	Undefined Antiviral Association
3	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 7 of 18

Plate	Well	Contents	Item Number	Primary Target*
3	B1	<i>Unused</i>		
3	B2	Isoborneol	23177	HSV-1 Protein Synthesis
3	B3	Isoxanthohumol	23233	Antiviral
3	B4	AG-555	23247	Mo-MuLV Reverse Transcriptase
3	B5	RN-18	23331	HIV-1 Vif
3	B6	5-(Hydroxymethyl)-2'-deoxyuridine	23381	HSV-1 Pyrimidine, 2'-deoxy Ribonucleoside Kinase
3	B7	20(S)-Ginsenoside Rh ₂	23390	P-glycoprotein, HIV
3	B8	Didanosine	23715	HIV Replication
3	B9	Hydroxyurea	23725	HIV-1 Replication
3	B10	Peramivir	23765	Influenza Neuraminidase
3	B11	Docosanol	23793	Broad-spectrum Antiviral (lipid-enveloped viruses)
3	B12	<i>Unused</i>		
3	C1	<i>Unused</i>		
3	C2	L-Valacyclovir (hydrochloride)	23801	Prodrug, Broad-spectrum Antiviral
3	C3	Famciclovir	23834	Prodrug, Herpes Viruses Replication
3	C4	Delavirdine (mesylate)	24026	HIV-1 Reverse Transcriptase
3	C5	Phenytoin	24037	Undefined Antiviral Association
3	C6	Hesperadin	24199	Influenza
3	C7	MST-312	24301	HSV replication
3	C8	Brequinar	24445	Undefined Antiviral Association
3	C9	Catalpol	24925	Undefined Antiviral Association
3	C10	Isoquercetin	24926	Influenza
3	C11	Aucubin	24928	Undefined Antiviral Association
3	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 8 of 18

Plate	Well	Contents	Item Number	Primary Target*
3	D1	<i>Unused</i>		
3	D2	Chicoric Acid	24960	HIV Integrase, HIV Entry
3	D3	3,5-Dicaffeoylquinic Acid	24964	HIV-1 Integrase
3	D4	Verbascoside	24965	Undefined Antiviral Association
3	D5	Senoside A	24969	Undefined Antiviral Association
3	D6	Saikosaponin A	25026	Undefined Antiviral Association
3	D7	Chelidonine	25097	Undefined Antiviral Association
3	D8	Theaflavin	25129	Undefined Antiviral Association
3	D9	Luteoloside	25141	Enterovirus 71
3	D10	Trilobatin	25837	HIV-1
3	D11	Imperatorin	26070	Antiviral
3	D12	<i>Unused</i>		
3	E1	<i>Unused</i>		
3	E2	GS-7340	26214	Prodrug, HIV-1 Replication
3	E3	Doravirine	26405	HIV Reverse Transcriptase
3	E4	Bergenin	26406	Undefined Antiviral Association
3	E5	Schisantherin B	26716	Undefined Antiviral Association
3	E6	(±)-BAY 41-4109	26907	HBV Replication
3	E7	Cabotegravir	27215	HIV-1 Integrase
3	E8	20(R)-Ginsenoside Rh ₂	27300	GHV Replication
3	E9	Oroxlylin A	27363	Influenza
3	E10	Des(benzylpyridyl) Atazanavir	27372	Metabolite of Atazanavir
3	E11	Scutellarin	27461	Undefined Antiviral Association
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 9 of 18

Plate	Well	Contents	Item Number	Primary Target*
3	F1	<i>Unused</i>		
3	F2	Schisandrin C	27475	Undefined Antiviral Association
3	F3	Pibrentasvir	27546	HCV Replication
3	F4	Columbianadin	27661	HIV-1 Replication
3	F5	Derenofylline	27666	Zika Virus Replication
3	F6	Narcissoside	27759	Epstein-Barr Virus Early Antigen
3	F7	Betulinic Aldehyde	27810	NS5 RNA-dependent RNA Polymerase, Dengue Virus Type 2 Replication
3	F8	Cridanimod (sodium salt)	27884	Broad-spectrum Antiviral
3	F9	4,5-Dicaffeoylquinic Acid	27892	HIV-1 Integrase, HIV-1 Replication
3	F10	Amenamevir	27921	HSV-1 Helicase-Primase Complex
3	F11	Glecaprevir	27934	HCV NS3/4A, HCV Replication
3	F12	<i>Unused</i>		
3	G1	<i>Unused</i>		
3	G2	Paritaprevir	28301	HCV NS3/4A, HCV Replication
3	G3	Adefovir dipivoxil	28373	Prodrug, HBV Replication
3	G4	PHA-680632	28381	Undefined Antiviral Association
3	G5	Coptisine (chloride)	28424	Undefined Antiviral Association
3	G6	ZK 756326 (hydrochloride)	28454	CCR8, HIV
3	G7	Pleconaril	28461	VP1 Capsid Protein, Rhinovirus, Enterovirus 71 Replication
3	G8	JCP251	28712	HIV Protease
3	G9	N ⁶ -Methyladenosine	28833	Undefined Antiviral Association
3	G10	RSV-604	29041	RSV Replication
3	G11	Fangchinoline	29243	HIV
3	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



Plate	Well	Contents	Item Number	Primary Target*
3	H1	<i>Unused</i>		
3	H2	Merimepodib	29351	Broad-spectrum Antiviral
3	H3	MK-8245	29421	HCV Replication
3	H4	Brivudine	29448	Herpes Viruses Replication
3	H5	Sophocarpine	29598	Undefined Antiviral Association
3	H6	Myriocin	63150	HCV Replication
3	H7	Miltefosine	63280	Undefined Antiviral Association
3	H7	U-0126	70970	Undefined Antiviral Association
3	H9	5-Aminolevulinic Acid (hydrochloride)	9001902	Undefined Antiviral Association
3	H10	Maslinic Acid	10009645	HIV-1 Replication
3	H11	SAHA	10009929	Undefined Antiviral Association
3	H12	<i>Unused</i>		
4	A1	<i>Unused</i>		
4	A2	Ursolic Acid	10072	Undefined Antiviral Association
4	A3	BSI-201	11304	PARP
4	A4	14-deoxy-11,12-didehydro Andrographolide	11670	Antiviral
4	A5	Asiatic Acid	11818	Undefined Antiviral Association
4	A6	18 β -Glycyrrhetic Acid	11845	Undefined Antiviral Association
4	A7	Gallic Acid	11846	Undefined Antiviral Association
4	A8	Trigonelline (chloride)	11904	Undefined Antiviral Association
4	A9	Punicalagin	13069	Broad-spectrum Antiviral
4	A10	D609 (potassium salt)	13307	Antiviral
4	A11	Arcyriaflavin A	13516	CMV Replication
4	A12	<i>Unused</i>		
4	B1	<i>Unused</i>		
4	B2	Tizoxanide	13693	Influenza, HCV, HBV
4	B3	Entecavir (hydrate)	13831	HBV, Hepadnaviruses
4	B4	Deferoxamine (mesylate)	14595	Undefined Antiviral Association
4	B5	Sophorose	14717	Virucidal
4	B6	<i>Unused</i>		
4	B7	Probucol	15043	Undefined Antiviral Association
4	B8	PF-429242 (hydrochloride)	15140	Dengue Virus
4	B9	PSI-7977	15402	HCV RNA Polymerase
4	B10	Indinavir (sulfate)	15150	HIV-1 Protease
4	B11	Galangin	15948	Undefined Antiviral Association
4	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 11 of 18

Plate	Well	Contents	Item Number	Primary Target*
4	C1	<i>Unused</i>		
4	C2	Zalcitabine	16019	Reverse Transcriptase
4	C3	Dapivirine	16247	HIV-1 Reverse Transcriptase
4	C4	7,8-Dihydroxyflavone	16946	Undefined Antiviral Association
4	C5	Chrysin	17402	Undefined Antiviral Association
4	C6	Hesperidin	18646	Undefined Antiviral Association
4	C7	Rutin (hydrate)	19868	Undefined Antiviral Association
4	C8	Methyl Gallate	19951	Undefined Antiviral Association
4	C9	Polydatin	21246	Undefined Antiviral Association
4	C10	Amantadine (hydrochloride)	21364	Influenza A M2 Proton Channel
4	C11	Timosaponin AIII	22236	Undefined Antiviral Association
4	C12	<i>Unused</i>		
4	D1	<i>Unused</i>		
4	D2	<i>Unused</i>		
4	D3	Mizoribine	23128	HCV RNA Replication
4	D4	T-705	23384	RNA-dependent RNA Polymerase, RNA Viruses
4	D5	Cobicistat	23433	CYP3A
4	D6	Rimantadine (hydrochloride)	23838	Matrix Protein 2
4	D7	Bitertanol	24054	Undefined Antiviral Association
4	D8	(-)-Galocatechin Gallate	24962	Undefined Antiviral Association
4	D9	3,4-Dicaffeoylquinic Acid	25053	Influenza
4	D10	Gentiopicroin	25102	Undefined Antiviral Association
4	D11	TAK-779	26744	HIV-1
4	D12	<i>Unused</i>		
4	E1	<i>Unused</i>		
4	E2	D-Pinitol	26815	Undefined Antiviral Association
4	E3	Geniposide	28281	Undefined Antiviral Association
4	E4	Cyclocytidine (hydrochloride)	29014	Undefined Antiviral Association
4	E5	20(S)-Protopanaxatriol	29073	Undefined Antiviral Association
4	E6	Cyclosporin C	29152	Vaccinia Virus Replication
4	E7	(±)-Catechin (hydrate)	29175	HCV RNA-dependent RNA Polymerase
4	E8	S-Acetyl-L-glutathione	29624	HSV-1 Replication
4	E9	N-Nonyldeoxyojirimycin	29698	Japanese Encephalitis Virus and Dengue Virus Type 2 Replication
4	E10	Geraniin	29924	HSV-1 and HSV-2 Replication
4	E11	OSW-1	30310	Corona Virus Replication
4	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 12 of 18

Plate	Well	Contents	Item Number	Primary Target*
4	F1	<i>Unused</i>		
4	F2	Remdesivir	30354	Prodrug, MERS-CoV and SARS-CoV
4	F3	GS-441524	30469	Active Metabolite, Broad-spectrum Antiviral
4	F4	Aspirin	70260	Undefined Antiviral Association
4	F5	tetramethyl Nordihydroguaiaretic Acid	70302	HSV Replication
4	F6	Baicalein	70610	Undefined Antiviral Association
4	F7	<i>trans</i> -Resveratrol	70675	Undefined Antiviral Association
4	F8	U-73122	70740	HIV-1 Integrase
4	F9	Chlorogenic Acid	70930	Influenza
4	F10	(-)-Epigallocatechin Gallate	70935	Undefined Antiviral Association
4	F11	Curcumin	81025	Undefined Antiviral Association
4	F12	<i>Unused</i>		
4	G1	<i>Unused</i>		
4	G2	Palmitoyl Ethanolamide	90350	Undefined Antiviral Association
4	G3	O-11	90510	HIV-1 Replication
4	G4	Pepstatin A	9000469	Influenza
4	G5	Resibufogenin	9001348	Undefined Antiviral Association
4	G6	MIV-150	9002360	Reverse Transcriptase
4	G7	KRIBB11	9002528	Dengue Virus
4	G8	Gemcitabine (hydrochloride)	9003096	Prodrug, Broad-spectrum Antiviral
4	G9	Luteolin	10004161	Undefined Antiviral Association
4	G10	Quercetin (hydrate)	10005169	Undefined Antiviral Association
4	G11	DL- α -Lipoic Acid	10005728	Undefined Antiviral Association
4	G12	<i>Unused</i>		
4	H1	<i>Unused</i>		
4	H2	Hesperetin	10006084	Undefined Antiviral Association
4	H3	CAY10567	10010233	HCV NS5B RNA-dependent RNA Polymerase
4	H4	Triciribine	10010237	Undefined Antiviral Association
4	H5	SB-415286	10010247	Undefined Antiviral Association
4	H6	AG-17	10010248	Undefined Antiviral Association
4	H7	Apigenin	10010275	Undefined Antiviral Association
4	H7	DRB	10010302	HIV Transcription
4	H9	Rolipram	10011132	Influenza
4	H10	NH125	10011250	Undefined Antiviral Association
4	H11	Plerixafor (hydrochloride hydrate)	10011332	HIV-1 Replication
4	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 13 of 18

Plate	Well	Contents	Item Number	Primary Target*
5	A1	<i>Unused</i>		
5	A2	GNF-2	16253	Undefined Antiviral Association
5	A3	GNF-5	16254	Undefined Antiviral Association
5	A4	SRPIN340	16284	HCV Replication
5	A5	PSI-6206	18391	Negative Control, HCV NS5B
5	A6	BMS 626529	18413	HIV-1 Attachment
5	A7	ABT-333	18482	RNA-directed RNA Polymerase
5	A8	(-)-Catechin	18644	Undefined Antiviral Association
5	A9	Ro 90-7501	18881	Vesicular Stomatitis Virus
5	A10	GS-9620	19628	TLR7, HIV-1, HBV
5	A11	Cepharanthine	19648	Antiviral
5	A12	<i>Unused</i>		
5	B1	<i>Unused</i>		
5	B2	Cinobufagin	19844	Undefined Antiviral Association
5	B3	Vacuolin-1	20425	Undefined Antiviral Association
5	B4	RO8191	20449	HCV Replication
5	B5	Foscarnet (sodium salt)	20774	DNA and RNA Polymerases, Reverse Transcriptase
5	B6	3',4',7-Trihydroxyisoflavone	20930	Undefined Antiviral Association
5	B7	PA-457	21339	HIV-1 Maturation
5	B8	Trifluorothymidine	21366	Thymidylate Synthase
5	B9	GSK2838232	21390	HIV-1 Maturation
5	B10	Gallotannin	21421	HCV Viral Entry
5	B11	SC-75741	21818	Influenza Replication
5	B12	<i>Unused</i>		
5	C1	<i>Unused</i>		
5	C2	Bithionol	21844	Host Caspases, Zika Virus
5	C3	Retro-2	21946	Retrograde Protein Trafficking, Antiviral
5	C4	Aloperine	22259	HIV-1 Replication and Envelope-mediated Cell-Cell Fusion
5	C5	G10	22353	Chikungunya Virus and Venezuelan Equine Encephalitis Virus Replication
5	C6	Usnic Acid	22585	Influenza
5	C7	Isoferulic Acid	22715	Undefined Antiviral Association
5	C8	Pirodavir	22922	Broad-spectrum Antipicornaviral
5	C9	2'-C-Methyladenosine	23062	HCV NS5B, HCV Replication
5	C10	Sappanone A	23205	Influenza Neuraminidase
5	C11	LDC-4297	23398	Broad-spectrum Viral Replication Inhibitor
5	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 14 of 18

Plate	Well	Contents	Item Number	Primary Target*
5	D1	<i>Unused</i>		
5	D2	Daclatasvir	23730	HCV NS5A
5	D3	Trimethobenzamide (hydrochloride)	24145	Undefined Antiviral Association
5	D4	20-hydroxy Arachidic Acid	24652	Rhinovirus Coat Protein
5	D5	Scutellarein	24977	Undefined Antiviral Association
5	D6	1,5-Dicaffeoylquinic Acid	25050	HIV-1 Integrase
5	D7	5-Caffeoylquinic Acid	25054	Undefined Antiviral Association
5	D8	Ginsenoside Ro	25137	Sendai Virus (Murine Respirovirus)
5	D9	Orientin	25143	Parainfluenza Type 3
5	D10	Chebulagic Acid	25203	Broad-spectrum Antiviral
5	D11	PSB-603	25637	Zika Virus
5	D12	<i>Unused</i>		
5	E1	<i>Unused</i>		
5	E2	Daphnoretin	26408	RSV Replication
5	E3	Bictegravir	26532	HIV-1 Integrase
5	E4	Lomibuvir	26630	HCV RNA-dependent RNA Polymerase, HCV Replication
5	E5	Emtricitabine S-oxide	26654	HIV-1 and HBV Reverse Transcriptases
5	E6	3-Deazaadenosine	9000785	Antiviral
5	E7	Saquinavir (mesylate)	9001893	HIV Protease
5	E8	Flavopiridol (hydrochloride)	10009197	CMV and HIV-1 Promoters, HIV-1 Replication
5	E9	PIK-93	10009212	Enterovirus RNA Replication
5	E10	Inotilone	10010089	Antiviral
5	E11	1-Deoxynojirimycin (hydrochloride)	10011718	HIV
5	E12	<i>Unused</i>		
5	F1	<i>Unused</i>		
5	F2	(-)-Levamisole (hydrochloride)	14874	Immunomodulator
5	F3	<i>Unused</i>		
5	F4	Losartan Carboxylic Acid	15957	AT ₁ Receptor
5	F5	Pyronaridine (phosphate)	16139	Antimalarial
5	F6	Pyrimethamine	16472	Antimalarial
5	F7	Hydroxychloroquine (sulfate)	17911	Antimalarial
5	F8	Primaquine (phosphate)	18223	Antimalarial
5	F9	GS-5885	18519	NS5A, HCV
5	F10	Ivermectin B _{1a}	18768	Antiparasitic
5	F11	Losartan Carboxaldehyde	18855	AT ₁ Receptor
5	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 15 of 18

Plate	Well	Contents	Item Number	Primary Target*
5	G1	<i>Unused</i>		
5	G2	Lumefantrine	20678	Antimalarial
5	G3	Piperaquine	20799	Antimalarial
5	G4	Sulfadoxin	21488	Antimalarial
5	G5	Mefloquine (hydrochloride)	23665	Antimalarial
5	G6	Quinine	23958	Antimalarial
5	G7	Losartan (potassium salt)	10006594	AT ₁ Receptor
5	G8	Artemether	11815	Antimalarial
5	G9	E-64d	13533	SARS-CoV, SARS-CoV-2
5	G10	PD 150606	13859	Undefined Antiviral Association
5	G11	Leupeptin (hemisulfate)	14026	Protease Inhibitor
5	G12	<i>Unused</i>		
5	H1	<i>Unused</i>		
5	H2	Doxycycline (hyclate)	14422	Antimalarial
5	H3	DBZ	14627	γ-Secretase
5	H4	Nafamostat (mesylate)	14837	Serine Protease Inhibitor
5	H5	Z-AEVD-FMK	14987	Caspase Inhibitor
5	H6	Chymostatin	15114	Protease Inhibitor
5	H7	Z-Asp-CH2-DCB	15143	Protease Inhibitor
5	H8	<i>Unused</i>		
5	H9	P005091	15224	Protease Inhibitor
5	H10	WP1130	15227	Protease Inhibitor
5	H11	UCF 101	15550	Protease Inhibitor
5	H12	<i>Unused</i>		
6	A1	<i>Unused</i>		
6	A2	Amodiaquine	15954	Antimalarial
6	A3	Camostat (mesylate)	16018	SARS-CoV, SARS-CoV-2
6	A4	Ximelagatran	16862	Thrombin Inhibitor
6	A5	Dabigatran etexilate	17131	Protease Inhibitor
6	A6	Dabigatran	17133	Thrombin Inhibitor
6	A7	Carfilzomib	17554	Proteasome Inhibitor
6	A8	ML-323	18010	Protease Inhibitor
6	A9	Pheleuin	19269	Undefined Antiviral Association
6	A10	Dihydroartemisinin	19846	Antimalarial Metabolite
6	A11	Quinidine	20356	Undefined Antiviral Association
6	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 16 of 18

Plate	Well	Contents	Item Number	Primary Target*
6	B1	<i>Unused</i>		
6	B2	HZ-1157	23436	HCV NS3/4A
6	B3	Gabexate (mesylate)	23767	Protease Inhibitor
6	B4	Danoprevir	24185	HCV NS3/4A
6	B5	CA-074	24679	Cathepsin B
6	B6	APC 366 (trifluoroacetate salt)	25990	Tryptase Inhibitor
6	B7	Antipain (hydrochloride)	26705	Protease Inhibitor
6	B8	VLX1570	29721	Ubiquitinase Inhibitor
6	B9	2-cyano-Pyrimidine	10007121	Cathepsin K
6	B10	E-64	10007963	Cystein Protease Inhibitor
6	B11	Bortezomib	10008822	Proteasome Inhibitor
6	B12	<i>Unused</i>		
6	C1	<i>Unused</i>		
6	C2	Dipyridamole	18189	SARS-CoV-2 Replication
6	C3	<i>Unused</i>		
6	C4	<i>Unused</i>		
6	C5	<i>Unused</i>		
6	C6	<i>Unused</i>		
6	C7	<i>Unused</i>		
6	C8	<i>Unused</i>		
6	C9	<i>Unused</i>		
6	C10	<i>Unused</i>		
6	C11	<i>Unused</i>		
6	C12	<i>Unused</i>		
6	D1	<i>Unused</i>		
6	D2	<i>Unused</i>		
6	D3	<i>Unused</i>		
6	D4	<i>Unused</i>		
6	D5	<i>Unused</i>		
6	D6	<i>Unused</i>		
6	D7	<i>Unused</i>		
6	D8	<i>Unused</i>		
6	D9	<i>Unused</i>		
6	D10	<i>Unused</i>		
6	D11	<i>Unused</i>		
6	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 17 of 18

Plate	Well	Contents	Item Number	Primary Target*
6	E1	Unused		
6	E2	Unused		
6	E3	Unused		
6	E4	Unused		
6	E5	Unused		
6	E6	Unused		
6	E7	Unused		
6	E8	Unused		
6	E9	Unused		
6	E10	Unused		
6	E11	Unused		
6	E12	Unused		
6	F1	Unused		
6	F2	Unused		
6	F3	Unused		
6	F4	Unused		
6	F5	Unused		
6	F6	Unused		
6	F7	Unused		
6	F8	Unused		
6	F9	Unused		
6	F10	Unused		
6	F11	Unused		
6	F12	Unused		
6	G1	Unused		
6	G2	Unused		
6	G3	Unused		
6	G4	Unused		
6	G5	Unused		
6	G6	Unused		
6	G7	Unused		
6	G8	Unused		
6	G9	Unused		
6	G10	Unused		
6	G11	Unused		
6	G12	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 18 of 18

Plate	Well	Contents	Item Number	Primary Target*
6	H1	<i>Unused</i>		
6	H2	<i>Unused</i>		
6	H3	<i>Unused</i>		
6	H4	<i>Unused</i>		
6	H5	<i>Unused</i>		
6	H6	<i>Unused</i>		
6	H7	<i>Unused</i>		
6	H8	<i>Unused</i>		
6	H9	<i>Unused</i>		
6	H10	<i>Unused</i>		
6	H11	<i>Unused</i>		
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