

PRODUCT INFORMATION



Off-Patent Drug Screening Library (96-Well)

Item No. 43505 • Batch No. 0809975

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

Page 1 of 22

Plate	Well	Contents	Item Number	Primary Target*
1	A1	Unused		
1	A2	PKC 412	10459	Multiple Kinases
1	A3	Capecitabine	10487	Thymidylate Synthase
1	A4	Candesartan Cilexetil	10489	AT ₁ Receptors
1	A5	ITF 2357	11045	Class I and II HDACs
1	A6	Nifedipine	11106	Ca _v 1 Channels
1	A7	Diphenhydramine (hydrochloride)	11158	Histamine H ₁ , Muscarinic M ₁₋₅ Receptors
1	A8	Venlafaxine (hydrochloride)	11567	DAT, NET, SERT
1	A9	Letrozole	11568	Aromatase
1	A10	Bexarotene	11571	RXR
1	A11	Eprosartan (mesylate)	11578	AT ₁ Receptors
1	A12	Unused		
1	B1	Unused		
1	B2	Glipizide	11579	K _{ATP} Channels
1	B3	Fluconazole	11594	Lanosterol 14- α -Demethylase
1	B4	Olmesartan Medoxomil	11614	AT ₁ Receptors
1	B5	Telmisartan	11615	AT ₁ Receptors
1	B6	Tamoxifen (citrate)	11629	Estrogen Receptors
1	B7	5-Fluorocytosine	11635	DNA Synthesis
1	B8	Amphotericin B	11636	Ergosterol Modulator
1	B9	Brinzolamide	11660	CAII, CAIV
1	B10	Gemcitabine	11690	DNA (Intercalating Agent)
1	B11	Lonafarnib	11746	FTase
1	B12	Unused		
1	C1	Unused		
1	C2	Olanzapine	11937	Various CNS Receptors
1	C3	Irbesartan	11952	AT ₁ Receptors
1	C4	Triptolide	11973	Various CNS Receptors
1	C5	Anastrozole	11987	Aromatase
1	C6	Olopatadine (hydrochloride)	11999	Histamine H ₁ Receptors
1	C7	Clozapine	12059	5-HT, Histamine, α_1 -Adrenergic Receptors
1	C8	1-hydroxy Vitamin D ₂	12077	Vitamin D Receptor
1	C9	Glimepiride	12090	K _{ATP} Channels
1	C10	Epirubicin (hydrochloride)	12091	DNA (Intercalating Agent)
1	C11	Tranilast	13044	
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, 08/22/2025

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 22

Plate	Well	Contents	Item Number	Primary Target*
1	D1	<i>Unused</i>		
1	D2	Oxaliplatin	13106	DNA (Cross-linking Agent)
1	D3	Carboplatin	13112	DNA (Cross-linking Agent)
1	D4	PA-824	13142	
1	D5	Sunitinib (malate)	13159	FLK1, FLT3, PDGFR β
1	D6	Gefitinib	13166	EGFR
1	D7	Agomelatine	13203	MT, 5-HT _{2B} , 5-HT _{2C} Receptors
1	D8	MS-275	13284	HDAC
1	D9	Itraconazole	13288	Lanosterol 14- α -Demethylase
1	D10	Piroxicam	13368	COX, NSAID
1	D11	Trandolapril	13522	ACE
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	Topiramate	13623	Multiple Targets
1	E3	Lansoprazole	13627	H ⁺ /K ⁺ ATPase
1	E4	Risperidone	13629	Multiple Targets
1	E5	<i>Unused</i>		
1	E6	Alfuzosin (hydrochloride)	13648	α_1 -Adrenergic Receptors
1	E7	Adapalene	13655	RAR
1	E8	Nitazoxanide	13692	Pyruvate:Ferredoxin Oxidoreductase
1	E9	Misoprostol	13820	EP ₂ , EP ₃ Receptors
1	E10	Entecavir (hydrate)	13831	HBV Reverse Transcriptase
1	E11	Cyclophosphamide (hydrate)	13849	DNA (Alkylating Agent)
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	Ganciclovir	13853	DNA Polymerase
1	F3	Timolol (maleate)	13974	β -Adrenergic Receptors
1	F4	Azathioprine	13987	Purine Biosynthesis
1	F5	Mycophenolate Mofetil	13988	IMPDH
1	F6	Sildenafil Citrate	14008	PDE5
1	F7	Prostaglandin E ₂	14010	EP Receptors
1	F8	Tadalafil	14024	PDE5
1	F9	Clofarabine	14125	Deoxycytidine Kinase, DNA Polymerase
1	F10	Febuxostat	14127	Xanthine Oxidase
1	F11	Topotecan (hydrochloride)	14129	Topoisomerase I
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 22

Plate	Well	Contents	Item Number	Primary Target*
1	G1	<i>Unused</i>		
1	G2	Entacapone	14153	Catechol O-methyltransferase
1	G3	Daunorubicin (hydrochloride)	14159	Topoisomerase
1	G4	Acyclovir	14160	DNA Polymerase
1	G5	Temozolomide	14163	DNA (Alkylating Agent)
1	G6	Idarubicin (hydrochloride)	14176	Topoisomerase II
1	G7	B-HT 920 (hydrochloride)	14177	Dopamine D ₂ Receptors
1	G8	Valsartan	14178	AT ₁ Receptors
1	G9	Irinotecan (hydrochloride hydrate)	14180	Topoisomerase I
1	G10	Amlexanox	14181	IKKε, TBK1
1	G11	Memantine (hydrochloride)	14184	NMDA Receptors
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	Bicalutamide	14250	Androgen Receptors
1	H3	Fludarabine Phosphate	14251	DNA Synthesis
1	H4	Flumazenil	14252	GABA _A Receptors
1	H5	Rivastigmine (tartrate)	14270	AChE, BChE
1	H6	Ciprofloxacin (hydrochloride hydrate)	14286	Topoisomerase II, Topoisomerase IV
1	H7	Verapamil (hydrochloride)	14288	L-type Calcium Channels
1	H8	Spectinomycin (hydrochloride hydrate)	14324	30S Ribosome, Protein Synthesis
1	H9	Tetracycline (hydrochloride)	14328	30S Ribosome, Protein Synthesis
1	H10	Efavirenz	14412	Reverse Transcriptase
1	H11	5-Fluorouracil	14416	Thymidylate Synthase
1	H12	<i>Unused</i>		
2	A1	<i>Unused</i>		
2	A2	Ampicillin (sodium salt)	14417	Penicillin-binding Proteins
2	A3	Fluoxetine (hydrochloride)	14418	SERT
2	A4	Rifampicin	14423	RNA Polymerase
2	A5	Darifenacin (hydrobromide)	14424	Muscarinic M ₁₋₅ Receptors
2	A6	Teniposide	14425	Topoisomerase II
2	A7	Nimodipine	14573	Ca _v 1.2 Channels
2	A8	Diazoxide	14576	SUR1/K _{ir} 6.2
2	A9	9-cis Retinoic Acid	14587	RAR, RXR
2	A10	Deferoxamine (mesylate)	14595	Iron Chelator
2	A11	Bromocriptine (mesylate)	14598	Various CNS Receptors
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 22

Plate	Well	Contents	Item Number	Primary Target*
2	B1	<i>Unused</i>		
2	B2	Sumatriptan (succinate)	14600	5-HT _{1B} , 5-HT _{1D} , 5-HT _{1F} Receptors
2	B3	Dorzolamide (hydrochloride)	14616	Carbonic Anhydrase
2	B4	Bumetanide	14630	NKCC1
2	B5	Homoharringtonine	14631	Human 80S Ribosome
2	B6	Lonidamine	14640	Hexokinase
2	B7	Maraviroc	14641	CCR5
2	B8	Vildagliptin	14705	DPP-4
2	B9	Vandetanib	14706	Multiple Kinases
2	B10	Dimethyl fumarate	14714	Nrf2
2	B11	Posaconazole	14737	Lanosterol 14- α -Demethylase
2	B12	<i>Unused</i>		
2	C1	<i>Unused</i>		
2	C2	Moxifloxacin (hydrochloride)	14830	DNA Gyrase
2	C3	Ibuprofen	14832	PDE4
2	C4	Gemfibrozil	14835	PPAR α , PPAR γ
2	C5	Nafamostat (mesylate)	14837	Serine Protease
2	C6	Sertraline (hydrochloride)	14839	SERT
2	C7	Mitoxantrone (hydrochloride)	14842	DNA (Intercalating Agent), Topoisomerase II, HIV Integrase
2	C8	Leflunomide	14860	AhR, DHODH
2	C9	Sildenafil	14866	α_{1A} -Adrenergic Receptors
2	C10	Aprepitant	14867	NK1
2	C11	Loperamide (hydrochloride)	14875	μ -Opioid Receptors
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	Pentostatin	14878	ADA
2	D3	Omeprazole	14880	H ⁺ /K ⁺ ATPase
2	D4	Amlodipine (besylate)	14886	L-type Calcium Channels
2	D5	Meloxicam	14906	COX-2, NSAID
2	D6	Finasteride	14938	5 α -Reductase Type II
2	D7	Rabeprazole (sodium salt)	14939	H ⁺ /K ⁺ ATPase
2	D8	<i>Unused</i>		
2	D9	Clenbuterol (hydrochloride)	14985	β_2 -Adrenergic Receptors
2	D10	Metyrapone	14994	CYP450
2	D11	Paroxetine (hydrochloride)	14998	SERT
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 22

Plate	Well	Contents	Item Number	Primary Target*
2	E1	<i>Unused</i>		
2	E2	Ondansetron (hydrochloride hydrate)	14999	5-HT ₃ Receptors
2	E3	Azithromycin	15004	Bacterial 50S Ribosomal Subunit
2	E4	Bazedoxifene (acetate)	15005	ER α , ER β
2	E5	Doxorubicin (hydrochloride)	15007	DNA (Intercalating Agent), Topoisomerase II
2	E6	Exemestane	15008	Aromatase
2	E7	(\pm)-lbutilide (hemifumarate)	15011	I _{Kr}
2	E8	Linezolid	15012	Bacterial 50S Ribosomal Subunit
2	E9	Methylprednisolone	15013	Glucocorticoid, Mineralocorticoid Receptors
2	E10	Miglitol	15014	α -Glucosidase
2	E11	Prazosin (hydrochloride)	15023	α_1 -Adrenergic, MT ₃ Receptors
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	Sulfasalazine	15025	
2	F3	Tigecycline	15026	Bacterial 30S Ribosomal Subunit
2	F4	Tolterodine (tartrate)	15027	Muscarinic M ₁₋₅ Receptors
2	F5	Varenicline (tartrate)	15030	nAChRs (α 7)
2	F6	Cilostazol	15035	PDE3A
2	F7	Reboxetine (mesylate)	15038	NET
2	F8	Dofetilide	15045	I _{Kr}
2	F9	Trifluoperazine (hydrochloride)	15068	α_{1A} -Adrenergic, Dopamine D ₂ Receptors
2	F10	Nevirapine	15117	Reverse Transcriptase
2	F11	<i>Unused</i>		
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	Roflumilast	15141	PDE4B, PDE4D
2	G3	Nelfinavir (mesylate)	15144	HIV-1 Protease
2	G4	Abiraterone Acetate	15148	CYP17A1
2	G5	Indinavir (sulfate)	15150	HIV-1 Protease
2	G6	Ketoconazole	15212	Lanosterol 14- α -Demethylase
2	G7	Amiodarone (hydrochloride)	15213	hERG
2	G8	Clotrimazole	15278	Lanosterol 14- α -Demethylase
2	G9	Minoxidil	15302	COX-1, K _{ATP} Channels
2	G10	Captopril	15313	ACE
2	G11	Auranofin	15316	Thioredoxin Reductase
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 22

Plate	Well	Contents	Item Number	Primary Target*
2	H1	<i>Unused</i>		
2	H2	Ebastine	15372	Histamine H ₁ Receptors
2	H3	Budesonide	15407	Glucocorticoid Receptors
2	H4	Carvedilol	15418	β-Adrenergic Receptors
2	H5	Miconazole (nitrate)	15420	Lanosterol 14-α-Demethylase
2	H6	Carbamazepine	15422	Voltage-gated Sodium Channels
2	H7	Ticagrelor	15425	P2Y ₁₂ Receptors
2	H8	Brimonidine (tartrate)	15426	α ₂ -Adrenergic Receptors
2	H9	Lamotrigine	15428	Na _v Channels
2	H10	Metoprolol (succinate)	15429	β ₁ -Adrenergic Receptors
2	H11	Idebenone	15475	Electron Transport Chain
2	H12	<i>Unused</i>		
3	A1	<i>Unused</i>		
3	A2	Oxaprozin	15476	COX-1, COX-2, NSAID
3	A3	Ramipril	15558	ACE
3	A4	Dexmedetomidine (hydrochloride)	15581	α ₂ -Adrenergic Receptors
3	A5	Dobutamine (hydrochloride)	15582	β-Adrenergic Receptors
3	A6	Naloxone (hydrochloride)	15594	Opioid Receptors
3	A7	Eplerenone	15616	Mineralocorticoid Receptors
3	A8	Fluvoxamine (maleate)	15617	Various CNS Receptors
3	A9	Loratadine	15625	Histamine H ₁ Receptors
3	A10	Voriconazole	15633	Lanosterol 14-α-Demethylase
3	A11	Ivabradine (hydrochloride)	15868	HCN Channels
3	A12	<i>Unused</i>		
3	B1	<i>Unused</i>		
3	B2	Moxalactam (sodium salt)	15869	Penicillin-binding Proteins
3	B3	Zolmitriptan	15871	5-HT _{1B} , 5-HT _{1D} Receptors
3	B4	Progesterone	15876	Progesterone Receptors
3	B5	Amitriptyline (hydrochloride)	15881	NET, SERT
3	B6	Cyclobenzaprine (hydrochloride)	15885	Muscarinic M ₁₋₃ , Histamine H ₁ , 5-HT ₂ Receptors
3	B7	Doxepin (hydrochloride)	15888	NET, SERT
3	B8	Maprotiline (hydrochloride)	15892	NET
3	B9	Nortriptyline (hydrochloride)	15904	NET, SERT
3	B10	Caspofungin (acetate)	15923	1,3-β-D-Glucan Synthase
3	B11	Edrophonium (chloride)	15928	AChE
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 22

Plate	Well	Contents	Item Number	Primary Target*
3	C1	<i>Unused</i>		
3	C2	<i>Unused</i>		
3	C3	Dutasteride	15956	5 α -Reductase
3	C4	Lofexidine (hydrochloride)	15989	α_2 -Adrenergic Receptors
3	C5	Cefoxitin (sodium salt)	15990	Penicillin-binding Proteins
3	C6	Salmeterol (xinafoate)	16009	β_2 -Adrenergic Receptors
3	C7	Camostat (mesylate)	16018	Serine Proteases
3	C8	Ciclopirox	16021	HIF-PH2
3	C9	Cefotaxime (sodium salt)	16040	Penicillin-binding Proteins, β -Lactamase
3	C10	Enalapril (maleate)	16041	ACE
3	C11	Rivaroxaban	16043	Factor Xa
3	C12	<i>Unused</i>		
3	D1	<i>Unused</i>		
3	D2	Meropenem (hydrate)	16068	Penicillin-binding Proteins
3	D3	Oseltamivir (phosphate)	16070	Neuraminidase
3	D4	Clomiphene (citrate)	16087	Estrogen Receptors
3	D5	Cefoperazone (sodium salt)	16113	Penicillin-binding Proteins
3	D6	Levosimendan	16128	Troponin C, K _{ATP} Channels, PDE3
3	D7	Chlorpromazine (hydrochloride)	16129	Various CNS Receptors
3	D8	Procarbazine (hydrochloride)	16133	DNA Methylation
3	D9	Ingenol-3-angelate	16207	PKC
3	D10	Phenoxybenzamine (hydrochloride)	16211	α -Adrenergic Receptors
3	D11	Benztropine (mesylate)	16214	Muscarinic M ₁ Receptors
3	D12	<i>Unused</i>		
3	E1	<i>Unused</i>		
3	E2	Ezetimibe	16331	NPC1L1
3	E3	Rifabutin	16468	Bacterial DNA-dependent RNA Polymerase
3	E4	Danazol	16471	Androgen Receptors
3	E5	Pyrimethamine	16472	DHFR
3	E6	Trimethoprim	16473	DHFR
3	E7	Rizatriptan (benzoate salt)	16475	5-HT _{1B} , 5-HT _{1D} Receptors
3	E8	Cyproterone Acetate	16622	Androgen Receptors
3	E9	Mozavaptan (hydrochloride)	16664	Vasopressin V ₂ Receptors
3	E10	Melphalan	16665	DNA (Alkylating Agent)
3	E11	Flupirtine (maleate)	16674	K _V 7
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 22

Plate	Well	Contents	Item Number	Primary Target*
3	F1	<i>Unused</i>		
3	F2	Unoprostone isopropyl ester	16681	FP Receptors
3	F3	Celiprolol (hydrochloride)	16715	β_1 -Adrenergic Receptors
3	F4	Palonosetron (hydrochloride)	16752	5-HT ₃ Receptors
3	F5	Deferasirox	16753	Iron Chelator
3	F6	Ribavirin	16757	Viral Polymerases
3	F7	PTC-124	16758	
3	F8	17-phenyl trinor Prostaglandin F _{2α} ethyl amide	16820	FP Receptors
3	F9	Ximelagatran	16862	Thrombin
3	F10	DL- α -Difluoromethylornithine (hydrochloride hydrate)	16889	Ornithine Decarboxylase
3	F11	Tenofovir Disoproxil (fumarate)	16922	HBV, HIV-1 Reverse Transcriptases
3	F12	<i>Unused</i>		
3	G1	<i>Unused</i>		
3	G2	Desloratadine	16931	Histamine H ₁ Receptors
3	G3	Ranitidine (hydrochloride)	16939	Histamine H ₂ Receptors
3	G4	Enoxacin	16956	DNA Gyrase, Topoisomerase IV
3	G5	Romidepsin	17130	Class I HDACs
3	G6	Moxidectin	17165	GABA _A Receptors
3	G7	Rebamipide	17186	Mucin
3	G8	Ifenprodil (hemitartrate)	17201	NMDA Receptors
3	G9	Esomeprazole (sodium salt)	17303	H ⁺ /K ⁺ ATPase
3	G10	Mirabegron	17319	β_3 -Adrenergic Receptors
3	G11	Solifenacin (succinate)	17320	Muscarinic M ₁₋₃ Receptors
3	G12	<i>Unused</i>		
3	H1	<i>Unused</i>		
3	H2	Temoporfin	17333	Pro-oxidant, Photosensitizer
3	H3	Verteporfin	17334	Endothelium, Photosensitizer
3	H4	Oxcarbazepine	17340	Calcium, Sodium, Potassium Channels
3	H5	Amoxapine	17347	NET, SERT, Various CNS Receptors
3	H6	Midodrine (hydrochloride)	17349	α_1 -Adrenergic Receptors
3	H7	Polmacoxib	17509	COX-2, CAI, CAII
3	H8	Felodipine	17535	Ca _v 1 Channels
3	H9	Nicardipine (hydrochloride)	17537	Ca _v 1 Channels
3	H10	Paricalcitol	17716	Vitamin D Receptor
3	H11	Sivelestat (sodium salt hydrate)	17779	Neutrophil Elastase
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 22

Plate	Well	Contents	Item Number	Primary Target*
4	A1	<i>Unused</i>		
4	A2	Stiripentol	17781	GABA _A Receptors
4	A3	Articaine (hydrochloride)	17883	Na _v Channels
4	A4	Pemiroloast (potassium salt)	17889	Histamine H ₁ Receptors
4	A5	Nitisinone	17924	4-Hydroxyphenylpyruvate Dioxygenase
4	A6	Dyphylline	17958	Adenosine A ₁₋₂ Receptors, PDE
4	A7	Felbamate	18000	GABA _A , NMDA Receptors
4	A8	Meisoindigo	18007	
4	A9	Triamcinolone acetonide	18026	Glucocorticoid Receptors
4	A10	Epinastine (hydrochloride)	18136	Histamine H ₁ Receptors
4	A11	Dipyridamole	18189	PDE5A
4	A12	<i>Unused</i>		
4	B1	<i>Unused</i>		
4	B2	Fexofenadine (hydrochloride)	18191	Histamine H ₁ Receptors
4	B3	Tolmetin (sodium salt hydrate)	18195	COX, NSAID
4	B4	Iloprost	18215	IP, EP ₁ Receptors
4	B5	Gestodene	18217	Progesterone Receptors
4	B6	Prostaglandin I ₂ (sodium salt)	18220	Platelets Aggregation
4	B7	Ciclesonide	18227	Glucocorticoid Receptors
4	B8	Beraprost (sodium salt)	18230	Platelets Aggregation
4	B9	(R)-Lansoprazole	18235	H ⁺ /K ⁺ ATPase
4	B10	Boceprevir	18379	HCV NS3/4A
4	B11	Azlocillin (sodium salt)	18424	Penicillin-binding Proteins
4	B12	<i>Unused</i>		
4	C1	<i>Unused</i>		
4	C2	Iopromide	18470	Contrast Agent
4	C3	Lamivudine	18514	HIV-1 Reverse Transcriptase
4	C4	Doxazosin (mesylate)	18633	α ₁ -Adrenergic Receptors
4	C5	Pentoxifylline	18720	Erythrocytes
4	C6	Oxiconazole (nitrate)	18724	Lanosterol 14-α-Demethylase
4	C7	Cimetidine	18743	Histamine H ₂ Receptors
4	C8	Cefonicid (sodium salt)	18745	Penicillin-binding Proteins
4	C9	Epothilone D	18765	Tubulin
4	C10	Rosuvastatin (calcium salt)	18813	HMG-CoA Reductase
4	C11	Nizatidine	18861	Histamine H ₂ 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 22

Plate	Well	Contents	Item Number	Primary Target*
4	D1	<i>Unused</i>		
4	D2	Ceftriaxone (sodium salt hydrate)	18866	Penicillin-binding Proteins
4	D3	Rufinamide	18870	CAVA, Na _v 1.1
4	D4	Mebendazole	18872	Tubulin
4	D5	Triclabendazole	18883	Helminths
4	D6	Droperidol	18884	Dopamine Receptors
4	D7	Methsuximide	19058	Calcium Channels
4	D8	Flavoxate (hydrochloride)	19096	Ca _v 1.2 Channels, Muscarinic Receptors
4	D9	Diflunisal	19187	COX, NSAID
4	D10	<i>Unused</i>		
4	D11	<i>Unused</i>		
4	D12	<i>Unused</i>		
4	E1	<i>Unused</i>		
4	E2	Flibanserin	19203	5-HT _{1A} , 5-HT _{2A} , Dopamine D ₄ Receptors
4	E3	Naftifine (hydrochloride)	19234	Squalene Epoxidase
4	E4	Fenofibric Acid	19262	PPARα
4	E5	Pitolisant	19307	Histamine H ₃ Receptors
4	E6	Repaglinide	19387	K _{ir} 6 Channels
4	E7	Clarithromycin	19455	23S rRNA
4	E8	Dirithromycin	19466	Bacterial 50S Ribosomal Subunit
4	E9	Cyproheptadine (hydrochloride hydrate)	19551	5-HT ₂ , Histamine H ₁ , Muscarinic Receptors
4	E10	Aliskiren (hemifumarate)	19640	Renin
4	E11	Bepidil (hydrochloride)	19645	Calcium, Sodium, Potassium Channels
4	E12	<i>Unused</i>		
4	F1	<i>Unused</i>		
4	F2	Cetirizine (hydrochloride)	19686	Histamine H ₁ Receptors
4	F3	Aztreonam	19784	Penicillin-binding Proteins
4	F4	Nalidixic Acid	19807	DNA Gyrase, Topoisomerase IV
4	F5	Etretinate	19878	
4	F6	Tolbutamide	19888	SUR1/K _{ir} 6.2
4	F7	Aripiprazole	19989	Dopamine D ₂ , -D _{2L} , -D ₃ , 5-HT _{1A} , 5-HT _{2B} Receptors
4	F8	Mirtazapine	19994	Various CNS Receptors
4	F9	Diltiazem (hydrochloride)	20079	Calcium Channels
4	F10	Lidocaine	20081	Ion Influx
4	F11	Econazole (nitrate)	20223	Lanosterol 14-α-Demethylase
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



Page 11 of 22

Plate	Well	Contents	Item Number	Primary Target*
4	G1	<i>Unused</i>		
4	G2	Rotigotine	20246	Dopamine D ₁₋₅ Receptors
4	G3	Nelarabine	20248	DNA Replication
4	G4	Nabumetone	20251	COX, NSAID
4	G5	Ketotifen (fumarate)	20303	Histamine H ₁ Receptors
4	G6	Rifapentine	20307	RNA Polymerase
4	G7	Triclosan	20342	FabI
4	G8	Flutamide	20359	Androgen Receptors
4	G9	Betamethasone	20363	Glucocorticoid Receptors
4	G10	<i>Unused</i>		
4	G11	Levofloxacin (hydrate)	20382	DNA Gyrase, Topoisomerase IV
4	G12	<i>Unused</i>		
4	H1	<i>Unused</i>		
4	H2	Deflazacort	20386	Glucocorticoid Receptors
4	H3	Flecainide (acetate)	20388	I _{Na} , I _{Kr}
4	H4	Ramelteon	20389	MT ₁ , MT ₂ Receptors
4	H5	Praziquantel	20508	Helminths
4	H6	Prednisone	20677	Glucocorticoid Receptors
4	H7	Acidinium (bromide)	20699	Muscarinic M ₁₋₅ Receptors
4	H8	Fluticasone Propionate	20703	Glucocorticoid Receptors
4	H9	Perphenazine	20735	Various CNS Receptors
4	H10	<i>Unused</i>		
4	H11	Prochlorperazine (maleate)	20742	Dopamine D ₂ Receptors
4	H12	<i>Unused</i>		
5	A1	<i>Unused</i>		
5	A2	Loxapine (succinate)	20760	Various CNS Receptors
5	A3	Piperacillin (sodium salt)	20766	Penicillin-binding Proteins
5	A4	Tolcapone	20768	Catechol O-methyltransferase
5	A5	<i>Unused</i>		
5	A6	Ticlopidine (hydrochloride)	20770	P2Y ₁₂ Receptors
5	A7	17 α -Estradiol	20776	Estrogen Receptors, 5 α -Reductase
5	A8	Etodolac	20833	COX, NSAID
5	A9	Asunaprevir	20835	HCV NS3/4A
5	A10	Acitretin	20853	CRABP-I, CRABP-II
5	A11	Toremifene (citrate)	20854	ER α , ER β
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



Page 12 of 22

Plate	Well	Contents	Item Number	Primary Target*
5	B1	<i>Unused</i>		
5	B2	Prednisolone	20866	Glucocorticoid Receptors
5	B3	Carbenicillin (sodium salt)	20871	Penicillin-binding Proteins
5	B4	Azelastine (hydrochloride)	20873	Histamine H ₁ Receptors
5	B5	Norethindrone	20941	Progesterone Receptors
5	B6	Etravirine	20946	Reverse Transcriptase
5	B7	Aminoglutethimide	20947	Aromatase
5	B8	Sulfamethazine	20976	Bacterial DHPS
5	B9	Mepyramine (maleate)	20978	Histamine H ₁ Receptors
5	B10	Brompheniramine (maleate)	20996	Histamine H ₁ Receptors
5	B11	Nisoldipine	20998	Calcium Channels
5	B12	<i>Unused</i>		
5	C1	<i>Unused</i>		
5	C2	(S)-(+)-Clopidogrel (sulfate)	21002	P2Y ₁₂ Receptors
5	C3	Levobunolol (hydrochloride)	21005	β-Adrenergic Receptors
5	C4	Nafcillin (sodium salt hydrate)	21008	Penicillin-binding Proteins
5	C5	Levocabastine (hydrochloride)	21014	Histamine H ₁ Receptors
5	C6	Tipranavir	21028	HIV Protease
5	C7	Etonogestrel	21062	Progesterone Receptors
5	C8	Metoprolol (tartrate)	21165	β ₁ -Adrenergic Receptors
5	C9	Bestatin	21217	Aminopeptidase
5	C10	Acetazolamide	21218	Carbonic Anhydrase
5	C11	Lomustine	21237	DNA (Alkylating Agent)
5	C12	<i>Unused</i>		
5	D1	<i>Unused</i>		
5	D2	Clobetasol Propionate	21251	Glucocorticoid Receptors
5	D3	Chlorpheniramine (maleate)	21253	Histamine H ₁ Receptors
5	D4	Tropium (chloride)	21279	Muscarinic M ₁₋₅ Receptors
5	D5	Alcaftadine	21290	Histamine H ₁ Receptors
5	D6	Terbutaline (hemisulfate)	21292	β ₂ -Adrenergic Receptors
5	D7	Indapamide	21308	K _v 1.1 Channels
5	D8	Torasemide	21312	Na ⁺ /Cl ⁻ Cotransporter
5	D9	Suprofen	21350	COX, NSAID
5	D10	<i>Unused</i>		
5	D11	Mometasone Furoate	21365	Glucocorticoid 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

Plate	Well	Contents	Item Number	Primary Target*
5	E1	<i>Unused</i>		
5	E2	Trifluorothymidine	21366	Thymidylate Synthase
5	E3	Rufloxacin (hydrochloride)	21413	DNA Gyrase
5	E4	(-)-Ropivacaine (hydrochloride hydrate)	21422	Sodium Channels
5	E5	Quinapril (hydrochloride)	21439	ACE
5	E6	Pindolol	21445	β_1 -, β_2 -Adrenergic Receptors
5	E7	Cefdinir	21521	Penicillin-binding Proteins
5	E8	Vilazodone (hydrochloride)	21547	5-HT _{1A} Receptors, SERT
5	E9	Megestrol Acetate	21749	Progesterone Receptors
5	E10	Tinidazole	21755	DNA (Oxidative Strand Breaks)
5	E11	Demeclocycline (hydrochloride)	21759	Bacterial 30S Ribosomal Subunit
5	E12	<i>Unused</i>		
5	F1	<i>Unused</i>		
5	F2	Nifurtimox	21784	
5	F3	Doxylamine (succinate)	21826	Histamine H ₁ Receptors
5	F4	<i>Unused</i>		
5	F5	Pemafibrate	22004	PPAR α
5	F6	Buspirone (hydrochloride)	22028	α -Adrenergic Receptors
5	F7	β -Estradiol 17-valerate	22221	Estrogen Receptors
5	F8	Sertaconazole (nitrate)	22232	Lanosterol 14- α -Demethylase
5	F9	Atomoxetine (hydrochloride)	22248	NET
5	F10	Pioglitazone (hydrochloride)	22263	PPAR γ
5	F11	Alosetron (hydrochloride)	22434	5-HT ₃ Receptors
5	F12	<i>Unused</i>		
5	G1	<i>Unused</i>		
5	G2	3-trifluoromethyl-N-Ethylamphetamine (hydrochloride)	22491	Various CNS Receptors
5	G3	Esmolol (hydrochloride)	22581	β_1 -Adrenergic Receptors
5	G4	Quetiapine (hemifumarate)	22603	Various CNS Receptors
5	G5	Hydrochlorothiazide	22671	SLC12A3
5	G6	Ofloxacin	22891	DNA Gyrase
5	G7	Guanfacine (hydrochloride)	22907	α_2 -Adrenergic Receptors
5	G8	<i>Unused</i>		
5	G9	Penciclovir	22918	DNA Polymerase
5	G10	Tiagabine (hydrochloride)	22926	GAT-1
5	G11	lloperidone	22957	Various CNS Receptors
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 22

Plate	Well	Contents	Item Number	Primary Target*
5	H1	<i>Unused</i>		
5	H2	Istradefylline	22958	Adenosine A _{2A} Receptors
5	H3	Cisatracurium (besylate)	22959	nAChRs
5	H4	Valrubicin	22964	DNA (Intercalating Agent), Topoisomerase II
5	H5	Malathion	22998	AChE
5	H6	Tetrahydrozoline (hydrochloride)	23004	α ₁ -Adrenergic Receptors
5	H7	Avanafil	23024	PDE5
5	H8	Clevidipine	23025	L-type Calcium Channels
5	H9	Luliconazole	23120	Lanosterol 14-α-Demethylase
5	H10	Hydralazine (hydrochloride)	23121	
5	H11	Pseudomonic Acid	23238	Bacterial Isoleucyl-tRNA Synthetase
5	H12	<i>Unused</i>		
6	A1	<i>Unused</i>		
6	A2	Citalopram (hydrobromide)	23252	SERT
6	A3	Clofazimine	23301	Mycobacteria
6	A4	Loteprednol etabonate	23305	Glucocorticoid Receptors
6	A5	Nateglinide	23320	K _{ATP} Channels
6	A6	Drospirenone	23347	Progesterone, Mineralocorticoid, Androgen Receptors
6	A7	(±)-Propranolol (hydrochloride)	23349	β-Adrenergic Receptors
6	A8	T-705	23384	RNA-dependent RNA Polymerase
6	A9	Tirofiban (hydrochloride hydrate)	23392	GPIIb/IIIa
6	A10	Methazolamide	23413	Carbonic Anhydrase
6	A11	Vesnarinone	23489	PDE3, hERG1
6	A12	<i>Unused</i>		
6	B1	<i>Unused</i>		
6	B2	Neostigmine (methyl sulfate)	23501	AChE
6	B3	Nitrofurantoin	23510	Bacterial Ribosomal Proteins
6	B4	Tasimelteon	23546	MT ₁ , MT ₂ Receptors
6	B5	Butoconazole (nitrate)	23554	Lanosterol 14-α-Demethylase
6	B6	Fluphenazine (hydrochloride)	23555	Dopamine D ₁ , D ₂ Receptors
6	B7	Tazarotene	23559	RAR
6	B8	Sulfamethoxazole	23613	DHPS
6	B9	Cefotetan (sodium salt)	23624	Penicillin-binding Proteins
6	B10	Isocarboxazid	23625	MAO
6	B11	Cefuroxime (sodium salt)	23627	Penicillin-binding Proteins
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 22

Plate	Well	Contents	Item Number	Primary Target*
6	C1	<i>Unused</i>		
6	C2	Fosfomycin Trometamol	23632	MurA
6	C3	Cefepime (hydrochloride hydrate)	23633	Penicillin-binding Proteins
6	C4	Mefenamic Acid	23650	COX-2, NSAID
6	C5	Ulipristal Acetate	23657	Progesterone Receptors
6	C6	Dichlorphenamide	23658	Carbonic Anhydrase
6	C7	Pralidoxime (chloride)	23666	AChE
6	C8	Ambrisentan	23669	ET _A Receptors
6	C9	6-Mercaptopurine (hydrate)	23675	DNA Synthesis
6	C10	Pralatrexate	23689	DHFR
6	C11	Besifloxacin (hydrochloride)	23692	Bacterial DNA Gyrase, Topoisomerase IV
6	C12	<i>Unused</i>		
6	D1	<i>Unused</i>		
6	D2	Rocuronium (bromide)	23698	nAChRs
6	D3	Albendazole	23705	Tubulin
6	D4	Mepenzolate (bromide)	23717	Muscarinic M ₂₋₃ Receptors
6	D5	Sulfadiazine	23719	DHPS
6	D6	Tiopronin	23720	Cystine
6	D7	Bepotastine (besylate)	23721	Histamine H ₁ Receptors
6	D8	Conivaptan (hydrochloride)	23728	Vasopressin V _{1A} , V ₂ Receptors
6	D9	Ixabepilone	23732	Tubulin
6	D10	Chlorambucil	23744	DNA (Alkylating Agent)
6	D11	Iohexol	23753	Contrast Agent
6	D12	<i>Unused</i>		
6	E1	<i>Unused</i>		
6	E2	Naratriptan (hydrochloride)	23754	5-HT _{1B} , 5-HT _{1D} Receptors
6	E3	Carteolol (hydrochloride)	23758	β-Adrenergic Receptors
6	E4	Dicloxacillin (sodium salt hydrate)	23770	Penicillin-binding Proteins
6	E5	Frovatriptan (succinate hydrate)	23771	5-HT _{1B} , 5-HT _{1D} Receptors
6	E6	Naphazoline (hydrochloride)	23774	α-Adrenergic Receptors
6	E7	Halobetasol Propionate	23781	Glucocorticoid Receptors
6	E8	Doxapram (hydrochloride hydrate)	23794	Carotid Chemoreceptors
6	E9	Butenafine (hydrochloride)	23797	Squalene Epoxidase
6	E10	Sulconazole (nitrate)	23800	Lanosterol 14-α-Demethylase
6	E11	L-Valacyclovir (hydrochloride)	23801	Viral DNA Polymerase
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 22

Plate	Well	Contents	Item Number	Primary Target*
6	F1	<i>Unused</i>		
6	F2	Atovaquone	23802	Mitochondrial Complex III
6	F3	Benazepril (hydrochloride)	23804	ACE
6	F4	Estropipate	23806	OATP1B1, Estrogen Receptors
6	F5	Oxybutynin (hydrochloride)	23825	Muscarinic Receptors
6	F6	Bisoprolol (hemifumarate)	23827	β_1 -Adrenergic Receptors
6	F7	Famotidine	23828	Histamine H ₂ Receptors
6	F8	Pyridostigmine (bromide)	23831	AChE, BChE
6	F9	Famciclovir	23834	DNA Polymerase
6	F10	Fenoprofen (calcium salt hydrate)	23835	NSAID
6	F11	Avobenzone	23836	UVA
6	F12	<i>Unused</i>		
6	G1	<i>Unused</i>		
6	G2	Milnacipran (hydrochloride)	23837	NET, SERT
6	G3	Rimantadine (hydrochloride)	23838	Influenza M2, HCV p7
6	G4	Efinaconazole	23839	Lanosterol 14- α -Demethylase
6	G5	Molindone (hydrochloride)	23841	Dopamine Receptors
6	G6	Dihydroergotamine (mesylate)	23847	5-HT _{1B} , 5-HT _{1D} Receptors
6	G7	Methocarbamol	23870	CAI
6	G8	Ropinirole (hydrochloride)	23871	Dopamine D ₂ Receptors
6	G9	Sertindole	23883	Various CNS Receptors
6	G10	Fosphenytoin (sodium salt)	23889	Na _v Channels
6	G11	Trimethadione	23893	
6	G12	<i>Unused</i>		
6	H1	<i>Unused</i>		
6	H2	<i>Unused</i>		
6	H3	Acetohexamide	23900	SUR1/K _{ir} 6.2
6	H4	Alclometasone Dipropionate	23901	Glucocorticoid Receptors
6	H5	Amcinonide	23903	Glucocorticoid Receptors
6	H6	Arformoterol (tartrate)	23935	β_2 -Adrenergic Receptors
6	H7	Benzonatate	23936	Na _v Channels
6	H8	Carbinoxamine (maleate)	23937	Histamine H ₁ Receptors
6	H9	Metaraminol (tartrate)	23938	α_1 -Adrenergic Receptors
6	H10	Cyclopentolate (hydrochloride)	23944	Muscarinic M ₁₋₃ Receptors
6	H11	Desoximetasone	23945	Glucocorticoid Receptors
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 22

Plate	Well	Contents	Item Number	Primary Target*
7	A1	<i>Unused</i>		
7	A2	Emedastine (fumarate)	23946	Histamine H ₁ Receptors
7	A3	Hexaminolevulinate (hydrochloride)	23949	Photosensitizer
7	A4	Levomilnacipran (hydrochloride)	23951	NET, SERT
7	A5	Nilutamide	23953	Androgen Receptors
7	A6	Medroxyprogesterone 17-acetate	23664	Progesterone Receptors
7	A7	Terconazole	23959	Lanosterol 14- α -Demethylase
7	A8	Cefamandole Nafate	23988	Penicillin-binding Proteins
7	A9	Isoflurane	23989	GABA _A Receptors
7	A10	(R)-Cetirizine (hydrochloride)	23992	Histamine H ₁ Receptors
7	A11	Sevoflurane	23996	GABA _A , Glycine Receptors
7	A12	<i>Unused</i>		
7	B1	<i>Unused</i>		
7	B2	Oxtriphylline	24006	Adenosine A ₁ , A ₂ Receptors, PDE
7	B3	Tamsulosin (hydrochloride)	24020	α_1 -Adrenergic Receptors
7	B4	Delavirdine (mesylate)	24026	HIV-1 Reverse Transcriptase
7	B5	Dipivefrin (hydrochloride)	24028	α_{1A} -, β_1 -, β_2 -Adrenergic Receptors
7	B6	Echothiophate (iodide)	24029	AChE
7	B7	Methylnaltrexone (bromide)	24072	μ -Opioid Receptors
7	B8	Zonisamide	24183	Sodium, Calcium Channels, Carbonic Anhydrase
7	B9	Sarpogrelate (hydrochloride)	24194	5-HT ₂ Receptors
7	B10	Prilocaine (hydrochloride)	24360	Na _v Channels
7	B11	<i>Unused</i>		
7	B12	<i>Unused</i>		
7	C1	<i>Unused</i>		
7	C2	Ceftibuten (hydrate)	25334	Penicillin-binding Proteins
7	C3	Betamethasone 17-valerate	25640	Glucocorticoid Receptors
7	C4	Chlormezanone	25716	
7	C5	Norfloxacin	25975	Bacterial DNA Gyrase, Topoisomerase IV
7	C6	Cefradine	26060	Penicillin-binding Proteins
7	C7	Cefalothin (sodium salt)	26063	Penicillin-binding Proteins
7	C8	Ticarcillin (sodium salt)	26065	Penicillin-binding Proteins
7	C9	Tioconazole	26069	Lanosterol 14- α -Demethylase
7	C10	Cilnidipine	26080	L-type, N-type Calcium Channels
7	C11	Ritodrine (hydrochloride)	26084	β_2 -Adrenergic Receptors
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 22

Plate	Well	Contents	Item Number	Primary Target*
7	D1	<i>Unused</i>		
7	D2	Pergolide (mesylate)	26085	Dopamine D ₁ , D ₂ Receptors
7	D3	Azilsartan	26091	AT ₁ Receptors
7	D4	Atenolol	26284	β ₁ -Adrenergic Receptors
7	D5	Betaxolol (hydrochloride)	26285	β ₁ -Adrenergic Receptors
7	D6	Labetalol (hydrochloride)	26289	α-, β-Adrenergic Receptors
7	D7	Furosemide	26298	NKCC1, NKCC2
7	D8	Ethacrynic Acid	26299	Na ⁺ /K ⁺ /2Cl ⁻ Symporter
7	D9	Cefmenoxime (hydrochloride)	26440	Penicillin-binding Proteins
7	D10	Lusutrombopag	26533	Thrombopoietin Receptor
7	D11	Scopolamine N-butyl (bromide)	27223	Muscarinic Receptors
7	D12	<i>Unused</i>		
7	E1	<i>Unused</i>		
7	E2	<i>Unused</i>		
7	E3	2',3',5'-Triacetyluridine	27445	Uridine, Dihydropyrimidine Dehydrogenase
7	E4	Trihexyphenidyl (hydrochloride)	27643	Muscarinic M ₁ Receptors
7	E5	Amenamivir	27921	HSV Helicase-Primase Complex
7	E6	Delafloxacin (meglumine)	27935	Bacterial DNA Gyrase, Topoisomerase IV
7	E7	<i>Unused</i>		
7	E8	Triamcinolone	28275	Glucocorticoid Receptors
7	E9	Adefovir dipivoxil	28373	DNA Synthesis
7	E10	Bilastine	28375	Histamine H ₁ Receptors
7	E11	Mavoxifafor	28439	CXCR4
7	E12	<i>Unused</i>		
7	F1	<i>Unused</i>		
7	F2	Palovarotene	28460	RARγ
7	F3	Pentagastrin	28546	CCK ₂ Receptor
7	F4	Benznidazole	29090	Protozoa
7	F5	Erdosteine	29177	
7	F6	Hydrocortisone Acetate	29390	Glucocorticoid Receptors
7	F7	Propentofylline	29431	PDE, Adenosine A ₁ , A ₂ Receptors, Adenosine Transporters
7	F8	<i>Unused</i>		
7	F9	Pirenzepine (hydrochloride)	29527	Muscarinic M ₁ Receptors
7	F10	<i>Unused</i>		
7	F11	Delamanid	29697	Mycolic Acid Synthesis
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 22

Plate	Well	Contents	Item Number	Primary Target*
7	G1	<i>Unused</i>		
7	G2	Fluticasone Furoate	29878	Glucocorticoid Receptors
7	G3	Isoconazole (nitrate)	30100	Heme Oxygenase
7	G4	Ilaprazole	30120	H ⁺ /K ⁺ ATPase
7	G5	Cefditoren Pivoxil	30315	Penicillin-binding Proteins
7	G6	Suplatast (tosylate)	30316	
7	G7	Udenafil	30373	PDE5
7	G8	Clevudine	30612	HBV DNA Polymerase
7	G9	Tebipenem Pivoxil	30613	Penicillin-binding Proteins
7	G10	<i>Unused</i>		
7	G11	Enoximone	30868	PDE3
7	G12	<i>Unused</i>		
7	H1	<i>Unused</i>		
7	H2	Estradiol Cypionate	31105	Estrogen Receptors
7	H3	Itopride (hydrochloride)	31183	AChE
7	H4	<i>Unused</i>		
7	H5	Prulifloxacin	31624	Bacterial DNA Gyrase, Topoisomerase IV
7	H6	Sparsentan	33429	AT ₁ , ET _A Receptors
7	H7	Gadoteridol	33577	Contrast Agent
7	H8	Furazolidone	34090	
7	H9	Picosulfate (sodium salt hydrate)	34566	
7	H10	Trabectedin	34662	DNA (Alkylating Agent), PXR
7	H11	<i>Unused</i>		
7	H12	<i>Unused</i>		
8	A1	<i>Unused</i>		
8	A2	Dabigatran etexilate (mesylate)	35248	Thrombin
8	A3	Odevixibat	35344	ASBT
8	A4	<i>Unused</i>		
8	A5	Talaporfin (sodium salt)	35722	Photosensitizer
8	A6	6 α -Methylprednisolone 21-hemisuccinate (sodium salt)	35742	Glucocorticoid Receptors
8	A7	6 α -Methylprednisolone 21-acetate	35799	Glucocorticoid Receptors
8	A8	Ceftizoxime (sodium salt)	36093	Penicillin-binding Proteins
8	A9	Anagliptin	36188	DPP-4
8	A10	Cefpirome (sulfate)	37398	Penicillin-binding Proteins
8	A11	(\pm)-Nefopam (hydrochloride)	38137	
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 22

Plate	Well	Contents	Item Number	Primary Target*
8	B1	<i>Unused</i>		
8	B2	Clonidine (hydrochloride)	38558	α_2 -Adrenergic Receptors
8	B3	Tizanidine (hydrochloride)	38559	α_2 -Adrenergic Receptors
8	B4	Beclomethasone Dipropionate	38880	Glucocorticoid Receptors
8	B5	D-Pantethine	39137	CoA
8	B6	Spirapril (hydrochloride)	39278	ACE
8	B7	Betamethasone Dipropionate	39499	Glucocorticoid Receptors
8	B8	<i>Unused</i>		
8	B9	Maribavir	39650	CMV pUL97
8	B10	Escitalopram (oxalate)	39841	SERT
8	B11	Fluphenazine decanoate	39903	Dopamine D ₁ , D ₂ Receptors
8	B12	<i>Unused</i>		
8	C1	<i>Unused</i>		
8	C2	Glycerol Phenylbutyrate	40832	
8	C3	Amibenonium (chloride)	40845	AChE
8	C4	Ozenoxacin	41837	Bacterial DNA Gyrase, Topoisomerase IV
8	C5	Triamcinolone hexacetonide	42600	Glucocorticoid Receptors
8	C6	<i>Unused</i>		
8	C7	Lornoxicam	70220	COX, NSAID
8	C8	(±)-Flurbiprofen	70250	COX, NSAID
8	C9	Indomethacin	70270	COX, NSAID
8	C10	(+)-Naproxen	70290	COX, NSAID
8	C11	Tolfenamic Acid	70480	COX-2, NSAID
8	C12	<i>Unused</i>		
8	D1	<i>Unused</i>		
8	D2	Diclofenac (sodium salt)	70680	COX, NSAID
8	D3	Ketorolac (tromethamine salt)	70690	COX, NSAID
8	D4	AM80	71770	RAR α
8	D5	Fluorometholone	9000157	Glucocorticoid Receptors
8	D6	Spironolactone	9000324	Aldosterone, Androgen, Mineralocorticoid Receptors
8	D7	Dronedarone (hydrochloride)	9000543	I _K
8	D8	<i>Unused</i>		
8	D9	Levetiracetam	9001820	SV2A
8	D10	Saquinavir (mesylate)	9001893	HIV Protease
8	D11	Seratrodast	9002014	TXA ₂ Receptor
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 22

Plate	Well	Contents	Item Number	Primary Target*
8	E1	<i>Unused</i>		
8	E2	Metronidazole	9002409	DNA Synthesis
8	E3	Dapoxetine (hydrochloride)	9002488	SERT
8	E4	Gemcitabine (hydrochloride)	9003096	DNA Synthesis
8	E5	(S)-Flurbiprofen	10004207	COX, NSAID
8	E6	Sulindac	10004386	COX, NSAID
8	E7	Fenofibrate	10005368	PPAR α
8	E8	Orlistat	10005426	Lipases
8	E9	Taflopust	10005440	FP Receptors
8	E10	Silybin	10006211	
8	E11	Fingolimod (hydrochloride)	10006292	S ₁ P Receptors
8	E12	<i>Unused</i>		
8	F1	<i>Unused</i>		
8	F2	17 β -Estradiol	10006315	ER α , ER β
8	F3	Mifepristone	10006317	Androgen, Glucocorticoid, Progesterone Receptors
8	F4	Levonorgestrel	10006318	Estrogen Receptors
8	F5	Norgestrel	10006319	Estrogen Receptors
8	F6	Quinestrol	10006320	Estrogen Receptors
8	F7	Ethynyl Estradiol	10006486	ER α , ER β
8	F8	Losartan (potassium salt)	10006594	AT ₁ Receptors
8	F9	(\pm)-Ketoprofen	10006661	COX, NSAID
8	F10	Zileuton	10006967	5-LO
8	F11	Zafirlukast	10008282	CysLT ₁
8	F12	<i>Unused</i>		
8	G1	<i>Unused</i>		
8	G2	Montelukast (sodium salt)	10008318	CysLT ₁
8	G3	Celecoxib	10008672	COX-2, NSAID
8	G4	Bortezomib	10008822	20S Proteasome
8	G5	Bezafibrate	10009145	PPAR
8	G6	SAHA	10009929	HDAC
8	G7	Lovastatin	10010338	HMG-CoA Reductase
8	G8	Pravastatin (sodium salt)	10010343	HMG-CoA Reductase
8	G9	Simvastatin	10010344	HMG-CoA Reductase
8	G10	Fulvestrant	10011269	ER α
8	G11	Terbinafine (hydrochloride)	10011619	Squalene Epoxidase
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

