

# PRODUCT INFORMATION



## Cellular Metabolism Screening Library (96-Well)

Item No. 33705 • Batch No. 0815640

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

Page 1 of 8

Plate	Well	Contents	Item Number	Primary Target*
1	A1	Unused		
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	$\Delta^6$ 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	Unused		
1	B1	Unused		
1	B2	Actinomycin D	11421	DNA Synthesis
1	B3	ABT-263	11500	Bcl-2, Bcl-xL, Bcl-W
1	B4	GSK1059615	11569	PI3K $\alpha$
1	B5	Phlorizin	11576	SGLT1, SGLT2
1	B6	Betulinic Acid	11686	TGR5
1	B7	Gemcitabine	11690	
1	B8	6-Gingerol	11707	TRPV1, TRPA1
1	B9	Artemisinin	11816	
1	B10	Artesunate	11817	GST
1	B11	6-Shogaol	11901	
1	B12	Unused		
1	C1	Unused		
1	C2	Dorsomorphin	11967	AMPK
1	C3	(+)-Etomoxir (sodium salt)	11969	CPT1
1	C4	Anastrozole	11987	CYP19A1
1	C5	Rottlerin	12006	Oxidative Phosphorylation
1	C6	CHIR99021	13122	GSK3
1	C7	JZL 184	13158	MAGL
1	C8	FK-866	13287	NAMPT
1	C9	Fatostatin (hydrobromide)	13562	SREBP
1	C10	YM-201636	13576	PIKfyve
1	C11	OU749	13804	GGT
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, 02/02/2026

#### 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 8

Plate	Well	Contents	Item Number	Primary Target*
1	D1	<i>Unused</i>		
1	D2	ML-265	13942	PKM2
1	D3	Methotrexate (hydrate)	13960	DHFR
1	D4	Rotenone	13995	Complex I
1	D5	AK-7	14004	SIRT2
1	D6	Febuxostat	14127	Xanthine Oxidase
1	D7	Fludarabine	14128	
1	D8	Plumbagin	14314	NOX
1	D9	AESBF (hydrochloride)	14321	NOX, Serine Proteases
1	D10	5-Fluorouracil	14416	Thymidylate Synthase
1	D11	Gossypol	14482	LDH, AKR1B1
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	Dorzolamide (hydrochloride)	14616	CAII, CAIV
1	E3	AGI-5198	14624	IDH1 R132H
1	E4	Bumetanide	14630	NKCC1, NKCC2
1	E5	AGI-6780	14639	IDH2 R140Q
1	E6	Lonidamine	14640	HK
1	E7	3-Nitropropionic Acid	14684	Complex II
1	E8	Dimethyl fumarate	14714	GSH
1	E9	TBCA	14831	CK2
1	E10	Pentostatin	14878	Adenosine Deaminase
1	E11	NB-598	14912	Squaline Epoxidase
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	Finasteride	14938	5 $\alpha$ -reductase II
1	F3	Metyrapone	14994	11 $\beta$ -Hydroxylase
1	F4	Linezolid	15012	Protein Synthesis
1	F5	Miglitol	15014	$\alpha$ -Glucosidase
1	F6	PF-04981517	15019	CYP3A4
1	F7	Torcetrapib	15028	CETP
1	F8	PF-429242 (hydrochloride)	15140	Site-1 Protease
1	F9	HQNO	15159	Complex III
1	F10	FCCP	15218	Oxidative Phosphorylation
1	F11	Atglistatin	15284	ATGL
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 8

Plate	Well	Contents	Item Number	Primary Target*
1	G1	<i>Unused</i>		
1	G2	Auranofin	15316	TxR1
1	G3	SF1670	15368	PTEN
1	G4	Ranolazine (hydrochloride)	15604	Sodium Channels
1	G5	CGP 37157	15611	Mitochondrial Sodium-Calcium Exchanger
1	G6	6-Thioguanine	15774	
1	G7	WWL113	16181	CES3, CES1F
1	G8	BMS 303141	16239	ACL
1	G9	PR-619	16276	DUBs
1	G10	TG101348	16289	JAK2, THTR2
1	G11	MitoTEMPO (hydrate)	16621	
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	U-104	16717	CAIX, CAXII
1	H3	Swainsonine	16860	Lysosomal $\alpha$ -Mannosidase, Golgi $\alpha$ -Mannosidase II
1	H4	DL- $\alpha$ -Difluoromethylornithine (hydrochloride hydrate)	16889	ODC
1	H5	UK 5099	16980	MPC
1	H6	CPI-613	16981	PDH, $\alpha$ -Ketoglutarate Dehydrogenase
1	H7	OGT 2115	17324	Heparanase
1	H8	Toyocamycin (hydrate)	17371	IRE1, PI3K, Rio1 Kinase
1	H9	GSK2656157	17372	PERK
1	H10	VPS34-IN-1	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, 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



Page 4 of 8

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	HIF-PH2
2	C3	WZB117	19900	Glut1
2	C4	AZD 3965	19912	MCT1
2	C5	GSK2194069	20022	FASN
2	C6	<i>Unused</i>		
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	Complex V
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	NCT-501	21162	ALDH1A1
2	D3	R-59-949	21221	DGK- $\alpha$
2	D4	Eliglustat (hemitartrate)	21487	GCS
2	D5	1-Deoxynojirimycin	21500	$\alpha$ -Glucosidase
2	D6	NLG-919 analog	21509	IDO, TDO
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	N/A
2	E3	6-Mercaptopurine (hydrate)	23675	Purine Synthesis
2	E4	ND-630	23961	ACC1, ACC2
2	E5	Orteronel	24191	CYP17A1
2	E6	ARN14988	24284	Acid Ceramidase
2	E7	Brequinar	24445	DHODH
2	E8	Raltitrexed	26079	Thymidylate Synthase
2	E9	CCF-642	27054	PDI
2	E10	Digitoxin	27825	HIF-1 $\alpha$
2	E11	DS44960156	28366	MTHFD2
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	iFSP1	29483	FSP1
2	F3	Liarozole	29491	CYP26A1
2	F4	Verinurad	29667	URAT1
2	F5	Alrestatin	29888	AKR1B1
2	F6	R162	30922	GDH1
2	F7	Pantothenate Kinase Inhibitor	31002	PanK
2	F8	Thioredoxin Reductase 1 Inhibitor 1	31307	TrxR1
2	F9	G6PDi-1	31484	G6PDH
2	F10	UGT8 Inhibitor 19	32723	UGT8
2	F11	Pachymic Acid	33180	HK2
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	N-Benzyl Linoleamide	33358	sEH
2	G3	LYS-006 (hydrochloride)	33706	LTA <sub>4</sub> hydrolase
2	G4	Enarodustat	33785	HIF-PH
2	G5	Meglutol	33832	HMG-CoA reductase
2	G6	DS18561882	33860	MTHFD2
2	G7	K6PC-5	34070	SPHK1
2	G8	LTI-291	34305	Glucocerebrosidase
2	G9	8-Azaadenosine	34622	ADAR1
2	G10	SH-42	34677	DHCR <sub>24</sub>
2	G11	JX06	34850	PDHK
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



Plate	Well	Contents	Item Number	Primary Target*
2	H1	<i>Unused</i>		
2	H2	Soticlestat	34928	Cholesterol 24-hydroxylase/CYP46A1
2	H3	BLT-1	34989	SCARB1
2	H4	JZP 430	35383	ABHD6
2	H5	SHIN1	35715	SHMT1, SHMT2
2	H6	Gemigliptin (tartrate)	36016	DPP-4
2	H7	Acetyl-CoA Synthase Inhibitor	36023	ACCS2
2	H8	JPH203 (hydrochloride)	36053	LAT1
2	H9	AG-270	36246	Methionine Adenosyltransferase 2A
2	H10	LY345899	36627	MTHFD
2	H11	GSK1940029	36704	Stearoyl-CoA Desaturase
2	H12	<i>Unused</i>		
3	A1	<i>Unused</i>		
3	A2	CTPI-2	36933	Mitochondrial Citrate Transporter/SLC25A1
3	A3	ARUK3001185	37113	NOTUM
3	A4	DUN41875	37151	Stearoyl-CoA Desaturase 1
3	A5	MDK8582	37307	sPLA <sub>2</sub>
3	A6	Defensamide	37318	SPHK1
3	A7	Verdiperstat	37333	MPO
3	A8	Imeglimin (hydrochloride)	37556	
3	A9	(R)-HTS-3	37563	LPCAT3
3	A10	FGTI-2734	37694	FTase, GGTase I
3	A11	Lipoxamycin (hemisulfate)	37747	Serine Palmitoyltransferase
3	A12	<i>Unused</i>		
3	B1	<i>Unused</i>		
3	B2	ML-198	37755	
3	B3	Dendrogenin A	37975	LXR, ChEH
3	B4	CAY10795	37981	15-PGDH
3	B5	Aryl piperazine 5k	38135	FATP1
3	B6	Prostaglandin E <sub>2</sub> Inhibitor 3	38316	mPGES-1
3	B7	AR-9281	38346	sEH
3	B8	UT-11	38580	mPGES-1
3	B9	ThioLox	38946	15-LO-1
3	B10	Ferroptocide	39005	Thioredoxin
3	B11	Oleoyl-L-carnitine (chloride)	39014	GlyT2
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

Plate	Well	Contents	Item Number	Primary Target*
3	C1	<i>Unused</i>		
3	C2	KYP-2047	39180	Prolyl Endopeptidase
3	C3	SLB1122168	39427	SPNS2
3	C4	VU534	39744	NAPE-PLD
3	C5	TP-0472993	39757	CYP4A11, CYP4F2
3	C6	Glutaminase C Inhibitor 11	39883	GAC
3	C7	<i>h</i> -e5'NT Inhibitor 5C	40266	Ecto-5'-nucleotidase
3	C8	ULK-101	40403	ULK1, ULK2
3	C9	Inaxaplin	40852	ApoL1
3	C10	K-604 (hydrochloride)	41119	ACAT-1
3	C11	BI-3231	41147	17 $\beta$ -HSD13
3	C12	<i>Unused</i>		
3	D1	<i>Unused</i>		
3	D2	BVT 14225	41224	11 $\beta$ -HSD1
3	D3	NNC-0640	41234	GCCR
3	D4	Glucagon Receptor Antagonist 2	41242	Glucagon Receptor
3	D5	BI-99179	41307	FAS-I
3	D6	SLU-PP-332	41719	ERR $\alpha$ , ERR $\beta$ , ERR $\gamma$
3	D7	Imirestat	41829	Aldose Reductase
3	D8	AZD 1656	41839	HK4
3	D9	BMS 466442	42075	Asc-1
3	D10	Insulin Sensitizer C59	42129	
3	D11	PSB-06126	42173	NTPDase 3
3	D12	<i>Unused</i>		
3	E1	<i>Unused</i>		
3	E2	UPGL00004	42323	GAC
3	E3	DMX-5084	42334	MAP4K4
3	E4	Talarozole	42372	CYP26
3	E5	TC-F 2	42395	FAAH
3	E6	Bz 423	42608	Complex V
3	E7	PLD <sub>1</sub> Inhibitor	42918	PLD <sub>1</sub>
3	E8	(E)-3-(2-Fluorophenyl)-N-(1H-indol-5-yl) acrylamide	43022	MAO-B, Nrf2
3	E9	IMT1	43029	POLRMT
3	E10	Z16078526	43072	UCP1
3	E11	2-Hydroxyisobutyric Acid	43901	
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



Plate	Well	Contents	Item Number	Primary Target*
3	F1	Unused		
3	F2	AZ-PFKFB3-67	43952	PFKFB3
3	F3	Semaglutide (trifluoroacetate salt)	44294	GLP-1R
3	F4	Myriocin	63150	Serine Palmitoyltransferase
3	F5	Miltefosine	63280	CTP:PCT
3	F6	SC-560	70340	COX-1
3	F7	Ebselen	70530	
3	F8	Chlorogenic Acid	70930	
3	F9	Rosiglitazone	71740	PPAR $\gamma$
3	F10	Diphenyleiiodonium (chloride)	81050	iNOS, eNOS
3	F11	NS 2028	81600	sGC
3	F12	Unused		
3	G1	Unused		
3	G2	Oleoyl Ethanolamide	90265	PPAR $\alpha$
3	G3	Vigabatrin	9000976	GABA-T
3	G4	STX-64	9003206	Steroid Sulfatase
3	G5	Orlistat	10005426	DAGL $\alpha$ , Lipases
3	G6	EUK 134	10006329	
3	G7	Ro 48-8071	10006415	OSC
3	G8	GW 0742	10006798	PPAR $\delta$
3	G9	AS-605240	10007707	PI3K $\alpha,\beta,\delta,\gamma$
3	G10	Celecoxib	10008672	COX-2
3	G11	Sphingosine Kinase Inhibitor 2	10009222	SPHK1
3	G12	Unused		
3	H1	Unused		
3	H2	6-Aminonicotinamide	10009315	G6PD
3	H3	( $\pm$ )-EX-527	10009798	SIRT1
3	H4	BMS309403	10010206	FABP4
3	H5	$\beta$ -cyano-L-Alanine	10010947	
3	H6	FTI-277 (hydrochloride)	10011667	FTase
3	H7	Eeyarestatin 1	10012609	ERAD
3	H8	A-922500	10012708	DGAT-1
3	H9	Unused		
3	H10	Unused		
3	H11	Unused		
3	H12	Unused		

\*Primary Target is based on available IC<sub>50</sub> and/or K<sub>i</sub> 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