

PRODUCT INFORMATION



Epigenetics Screening Library (96-Well)

Item No. 11076 • Batch No. 0605203

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

Page 1 of 6

Plate	Well	Contents	Item Number
1	A1	<i>Unused</i>	
1	A2	Sulforaphane	10496
1	A3	UNC669	10875
1	A4	SB939	10443
1	A5	PCI 34051	10444
1	A6	L- α -Hydroxyglutaric Acid (sodium salt)	11876
1	A7	Sirtinol	10523
1	A8	5-Nitroso-8-quinolinol	13856
1	A9	C646	10549
1	A10	JNJ-26481585	14088
1	A11	Garcinol	10566
1	A12	<i>Unused</i>	
1	B1	<i>Unused</i>	
1	B2	Plumbagin	14314
1	B3	Scriptaid	10572
1	B4	Suberohydroxamic Acid	10574
1	B5	Apicidin	10575
1	B6	LMK 235	14969
1	B7	UNC0321 (trifluoroacetate salt)	10582
1	B8	HDAC6 Inhibitor	15200
1	B9	Cl-Amidine (hydrochloride)	10599
1	B10	SP2509	15487
1	B11	JGB1741	10641
1	B12	<i>Unused</i>	
1	C1	<i>Unused</i>	
1	C2	A-366	16081
1	C3	I-BET762	10676
1	C4	UNC0638	10734
1	C5	OICR-9429	16095
1	C6	EPZ004777 (formate)	16173
1	C7	EPZ6438	16174
1	C8	PBIT	16272
1	C9	Zebularine	10975
1	C10	Delphinidin (chloride)	11012
1	C11	ITF 2357	11045
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, 05/27/2021

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 6

Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	Octyl- α -hydroxyglutarate	16365
1	D3	UNC0379	16400
1	D4	Methylstat (hydrate)	11091
1	D5	BVT 948	16615
1	D6	BG45	16918
1	D7	UMB-32	17123
1	D8	2,4-Pyridinedicarboxylic Acid (hydrate)	11138
1	D9	PFI-1	11155
1	D10	5-Azacytidine	11164
1	D11	SGI-1027	11165
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	Decitabine	11166
1	E3	I-BET151	11181
1	E4	(+)-JQ1	11187
1	E5	EPZ015666	17285
1	E6	BI-2536	17385
1	E7	BAZ2-ICR	17448
1	E8	Sodium 4-Phenylbutyrate	11323
1	E9	OG-L002	17471
1	E10	IOX1	11572
1	E11	MI-2 (hydrochloride)	11620
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	ML-324	17472
1	F3	Gemcitabine	11690
1	F4	GSK484 (hydrochloride)	17488
1	F5	Resminostat (hydrochloride)	17553
1	F6	NI-57	17662
1	F7	Daminozide	12033
1	F8	GSK-J1 (sodium salt)	12054
1	F9	PFI-4	17663
1	F10	GSK-J4 (hydrochloride)	12073
1	F11	Tasquinimod	17692
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
1	G1	<i>Unused</i>	
1	G2	CI-994	12084
1	G3	CPTH2 (hydrochloride)	12086
1	G4	MM-102	17699
1	G5	TC-E 5003	17718
1	G6	Butyrolactone 3	12095
1	G7	L002	17778
1	G8	Tenovin-1	13085
1	G9	Tenovin-6	13086
1	G10	BI-9564	17897
1	G11	RN-1 (hydrochloride)	18124
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	Anacardic Acid	13144
1	H3	AGK2	13145
1	H4	CAY10603	13146
1	H5	Splitomicin	13168
1	H6	CBHA	13172
1	H7	Mocetinostat	18287
1	H8	Oxamflatin	13176
1	H9	Salermide	13178
1	H10	A-196	18317
1	H11	MS049 (hydrochloride)	18348
1	H12	<i>Unused</i>	
2	A1	<i>Unused</i>	
2	A2	GSK591	18354
2	A3	Panobinostat	13280
2	A4	MS-275	13284
2	A5	MS023 (hydrochloride)	18361
2	A6	RG108	13302
2	A7	2',3',5'-triacetyl-5-Azacytidine	13373
2	A8	PRT4165	19093
2	A9	UNC0224	13631
2	A10	ORY-1001	19136
2	A11	EPZ020411	19160
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 6

Plate	Well	Contents	Item Number
2	B1	<i>Unused</i>	
2	B2	3-Deazaneplanocin A	13828
2	B3	Sinefungin	13829
2	B4	Pyroxamide	13870
2	B5	N-Oxalylglycine	13944
2	B6	GSK2879552	19403
2	B7	AMI-1 (sodium salt)	13965
2	B8	EPZ005687	13966
2	B9	SGC0946	13967
2	B10	UNC1215	13968
2	B11	Cyproheptadine (hydrochloride hydrate)	19551
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	GSK343	14094
2	C3	Bromosporine	14119
2	C4	GSK2801	14120
2	C5	SIRT1/2 Inhibitor IV	14407
2	C6	I-CBP112 (hydrochloride)	14468
2	C7	SGC-CBP30	14469
2	C8	UNC0642	14604
2	C9	AZD 5153	20864
2	C10	(R)-PFI-2 (hydrochloride)	14678
2	C11	HPOB	15066
2	C12	<i>Unused</i>	
2	D1	<i>Unused</i>	
2	D2	CPTH6 (hydrobromide)	19828
2	D3	PFI-3	15267
2	D4	JIB-04	15338
2	D5	CAY10683	15403
2	D6	GSK126	15415
2	D7	CPI-203	15479
2	D8	6-Thioguanine	15774
2	D9	HAT Inhibitor II	19835
2	D10	CeMMEC1	19956
2	D11	OTX015	15947
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 6

Plate	Well	Contents	Item Number
2	E1	<i>Unused</i>	
2	E2	PCI 24781	20059
2	E3	BML-278	20209
2	E4	CeMMEC13	20224
2	E5	EPZ5676	16175
2	E6	MC 1568	16265
2	E7	4'-bromo-Resveratrol	20541
2	E8	α -Hydroxyglutaric Acid (sodium salt)	16374
2	E9	S-(5'-Adenosyl)-L-methionine (tosylate)	16376
2	E10	RVX-208	16424
2	E11	Todalazine (hydrochloride)	20893
2	E12	<i>Unused</i>	
2	F1	<i>Unused</i>	
2	F2	LAQ824	16427
2	F3	HDAC3 Inhibitor	21057
2	F4	GSK-LSD1 (hydrochloride)	16439
2	F5	CXD101	21159
2	F6	RGFP966	16917
2	F7	BRD73954	16919
2	F8	NCH-51	21234
2	F9	UF010	21273
2	F10	RGFP109	21568
2	F11	Trichostatin A	89730
2	F12	<i>Unused</i>	
2	G1	<i>Unused</i>	
2	G2	CAY10398	89740
2	G3	RSC-133	9001839
2	G4	BML-210	10005019
2	G5	C-7280948	21601
2	G6	CAY10591	10009797
2	G7	EED226	22031
2	G8	SAHA	10009929
2	G9	Tranylcypromine (hydrochloride)	10010494
2	G10	CAY10722	22404
2	G11	PAOA	22942
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

