

PRODUCT INFORMATION



Epigenetics Screening Library (96-Well)

Item No. 11076 • Batch No. 0697870

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

Page 1 of 8

Plate	Well	Contents	Item Number
1	A1	<i>Unused</i>	
1	A2	SB939	10443
1	A3	PCI 34051	10444
1	A4	4-iodo-SAHA	10495
1	A5	Sulforaphane	10496
1	A6	Sirtinol	10523
1	A7	C646	10549
1	A8	Garcinol	10566
1	A9	Ellagic Acid	10569
1	A10	Scriptaid	10572
1	A11	Suberohydroxamic Acid	10574
1	A12	<i>Unused</i>	
1	B1	<i>Unused</i>	
1	B2	Apicidin	10575
1	B3	Cl-Amidine (hydrochloride)	10599
1	B4	F-Amidine (trifluoroacetate salt)	10610
1	B5	I-BET762	10676
1	B6	UNC0638	10734
1	B7	UNC669	10875
1	B8	CAY10669	10974
1	B9	Zebularine	10975
1	B10	Delphinidin (chloride)	11012
1	B11	ITF 2357	11045
1	B12	<i>Unused</i>	
1	C1	<i>Unused</i>	
1	C2	UNC0631	11084
1	C3	UNC0646	11085
1	C4	Methylstat (hydrate)	11091
1	C5	2,4-Pyridinedicarboxylic Acid (hydrate)	11138
1	C6	PFI-1	11155
1	C7	5-Azacytidine	11164
1	C8	SGI-1027	11165
1	C9	Decitabine	11166
1	C10	I-BET151	11181
1	C11	(+)-JQ1	11187
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, 01/5/2024

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



Page 2 of 8

Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	Sodium 4-Phenylbutyrate	11323
1	D3	IOX1	11572
1	D4	Tofacitinib (citrate)	11598
1	D5	MI-2 (hydrochloride)	11620
1	D6	Gemcitabine	11690
1	D7	Daminozide	12033
1	D8	GSK-J1 (sodium salt)	12054
1	D9	GSK-J4 (hydrochloride)	12073
1	D10	CI-994	12084
1	D11	CPTH2 (hydrochloride)	12086
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	Butyrolactone 3	12095
1	E3	Valproic Acid (sodium salt)	13033
1	E4	Tenovin-1	13085
1	E5	Tenovin-6 (hydrochloride)	13086
1	E6	<i>Unused</i>	
1	E7	Anacardic Acid	13144
1	E8	AGK2	13145
1	E9	CAY10603	13146
1	E10	Splitomicin	13168
1	E11	CBHA	13172
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	M 344	13174
1	F3	Oxamflatin	13176
1	F4	Salermide	13178
1	F5	Pimelic Diphenylamide 106	13212
1	F6	Panobinostat	13280
1	F7	MS-275	13284
1	F8	HNHA	13295
1	F9	RG108	13302
1	F10	2',3',5'-triacetyl-5-Azacytidine	13373
1	F11	UNC0224	13631
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
1	G1	<i>Unused</i>	
1	G2	Chidamide	13686
1	G3	3-Deazaneplanocin A	13828
1	G4	5-Nitroso-8-quinolinol	13856
1	G5	Pyroxamide	13870
1	G6	N-Oxalylglycine	13944
1	G7	WDR5-0103	13945
1	G8	AMI-1 (sodium salt)	13965
1	G9	EPZ005687	13966
1	G10	<i>Unused</i>	
1	G11	UNC1215	13968
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	AK-7	14004
1	H3	JNJ-26481585 (hydrochloride)	14088
1	H4	GSK343	14094
1	H5	Bromosporine	14119
1	H6	GSK2801	14120
1	H7	Plumbagin	14314
1	H8	SIRT1/2 Inhibitor IV	14407
1	H9	I-CBP112 (hydrochloride)	14468
1	H10	SGC-CBP30	14469
1	H11	UNC0642	14604
1	H12	<i>Unused</i>	
2	A1	<i>Unused</i>	
2	A2	UNC1999	14621
2	A3	(R)-PFI-2 (hydrochloride)	14678
2	A4	LMK 235	14969
2	A5	HPOB	15066
2	A6	HDAC6 Inhibitor	15200
2	A7	2-hexyl-4-Pentynoic Acid	15205
2	A8	PFI-3	15267
2	A9	JIB-04	15338
2	A10	CAY10683	15403
2	A11	GSK126	15415
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
2	B1	<i>Unused</i>	
2	B2	CPI-203	15479
2	B3	SP-2509	15487
2	B4	6-Thioguanine	15774
2	B5	Tubastatin A	15785
2	B6	OTX015	15947
2	B7	A-366	16081
2	B8	OICR-9429	16095
2	B9	EPZ004777 (formate)	16173
2	B10	EPZ6438	16174
2	B11	EPZ5676	16175
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	MC 1568	16265
2	C3	PBIT	16272
2	C4	Octyl- α -hydroxyglutarate	16365
2	C5	α -Hydroxyglutaric Acid (sodium salt)	16374
2	C6	UNC0379	16400
2	C7	RVX-208	16424
2	C8	CUDC-101	16426
2	C9	<i>Unused</i>	
2	C10	GSK-LSD1 (hydrochloride)	16439
2	C11	BVT 948	16615
2	C12	<i>Unused</i>	
2	D1	<i>Unused</i>	
2	D2	RGFP966	16917
2	D3	BG45	16918
2	D4	BRD73954	16919
2	D5	UMB-32	17123
2	D6	EPZ015666	17285
2	D7	BI-2536	17385
2	D8	BAZ2-ICR	17448
2	D9	OG-L002	17471
2	D10	ML-324	17472
2	D11	GSK484 (hydrochloride)	17488
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 8

Plate	Well	Contents	Item Number
2	E1	<i>Unused</i>	
2	E2	Resminostat (hydrochloride)	17553
2	E3	NI-57	17662
2	E4	PFI-4	17663
2	E5	Tasquinimod	17692
2	E6	MM-102	17699
2	E7	TC-E 5003	17718
2	E8	Sephin1	17757
2	E9	L002	17778
2	E10	BI-9564	17897
2	E11	RN-1 (hydrochloride)	18124
2	E12	<i>Unused</i>	
2	F1	<i>Unused</i>	
2	F2	Mocetinostat	18287
2	F3	MCB-613	18315
2	F4	A-196	18317
2	F5	MS049 (hydrochloride)	18348
2	F6	GSK591	18354
2	F7	MS023 (hydrochloride)	18361
2	F8	PRT4165	19093
2	F9	EPZ020411	19160
2	F10	GSK2879552	19403
2	F11	Cyproheptadine (hydrochloride hydrate)	19551
2	F12	<i>Unused</i>	
2	G1	<i>Unused</i>	
2	G2	CPTH6 (hydrobromide)	19828
2	G3	HAT Inhibitor II	19835
2	G4	CeMMEC1	19956
2	G5	PCI 24781	20059
2	G6	BML-278	20209
2	G7	CeMMEC13	20224
2	G8	4'-bromo-Resveratrol	20541
2	G9	AZD 5153	20864
2	G10	Todralazine (hydrochloride)	20893
2	G11	HDAC3 Inhibitor	21057
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 8

Plate	Well	Contents	Item Number
2	H1	<i>Unused</i>	
2	H2	NCH-51	21234
2	H3	UF010	21273
2	H4	RGFP109	21568
2	H5	C-7280948	21601
2	H6	EED226	22031
2	H7	CAY10722	22404
2	H8	PAOA	22942
2	H9	TMP-195	23242
2	H10	CM-272	25948
2	H11	Bufexamac	26068
2	H12	<i>Unused</i>	
3	A1	<i>Unused</i>	
3	A2	DS-437	29476
3	A3	NR-160	31792
3	A4	BMS-P5	33581
3	A5	Belinostat	34084
3	A6	Trotabresib	36292
3	A7	FB23-2	36606
3	A8	ML-792	36820
3	A9	FL-411	37547
3	A10	UNC8153 (trifluoroacetate salt)	38932
3	A11	Trichostatin A	89730
3	A12	<i>Unused</i>	
3	B1	<i>Unused</i>	
3	B2	CAY10398	89740
3	B3	RSC-133	9001839
3	B4	BML-210	10005019
3	B5	CAY10591	10009797
3	B6	(±)-EX-527	10009798
3	B7	SAHA	10009929
3	B8	Tranilcypromine (hydrochloride)	10010494
3	B9	BIX01294 (hydrochloride hydrate)	13124
3	B10	<i>Unused</i>	
3	B11	<i>Unused</i>	
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 8

Plate	Well	Contents	Item Number
3	C1	Unused	
3	C2	Unused	
3	C3	Unused	
3	C4	Unused	
3	C5	Unused	
3	C6	Unused	
3	C7	Unused	
3	C8	Unused	
3	C9	Unused	
3	C10	Unused	
3	C11	Unused	
3	C12	Unused	
3	D1	Unused	
3	D2	Unused	
3	D3	Unused	
3	D4	Unused	
3	D5	Unused	
3	D6	Unused	
3	D7	Unused	
3	D8	Unused	
3	D9	Unused	
3	D10	Unused	
3	D11	Unused	
3	D12	Unused	
3	E1	Unused	
3	E2	Unused	
3	E3	Unused	
3	E4	Unused	
3	E5	Unused	
3	E6	Unused	
3	E7	Unused	
3	E8	Unused	
3	E9	Unused	
3	E10	Unused	
3	E11	Unused	
3	E12	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 8 of 8

Plate	Well	Contents	Item Number
3	F1	Unused	
3	F2	Unused	
3	F3	Unused	
3	F4	Unused	
3	F5	Unused	
3	F6	Unused	
3	F7	Unused	
3	F8	Unused	
3	F9	Unused	
3	F10	Unused	
3	F11	Unused	
3	F12	Unused	
3	G1	Unused	
3	G2	Unused	
3	G3	Unused	
3	G4	Unused	
3	G5	Unused	
3	G6	Unused	
3	G7	Unused	
3	G8	Unused	
3	G9	Unused	
3	G10	Unused	
3	G11	Unused	
3	G12	Unused	
3	H1	Unused	
3	H2	Unused	
3	H3	Unused	
3	H4	Unused	
3	H5	Unused	
3	H6	Unused	
3	H7	Unused	
3	H8	Unused	
3	H9	Unused	
3	H10	Unused	
3	H11	Unused	
3	H12	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