

# PRODUCT INFORMATION



## Epigenetics Screening Library (96-Well)

Item No. 11076 • Batch No. 0522205

The composition of the library will always vary somewhat depending upon our inventory.

Page 1 of 5

Plate	Well	Contents	Item Number
1	A1	<i>Unused</i>	
1	A2	(+)-Abscisic Acid	10073
1	A3	3-amino Benzamide	10397
1	A4	SB939	10443
1	A5	PCI 34051	10444
1	A6	4-iodo-SAHA	10495
1	A7	Sirtinol	10523
1	A8	C646	10549
1	A9	Tubastatin A (trifluoroacetate salt)	10559
1	A10	Garcinol	10566
1	A11	Ellagic Acid	10569
1	A12	<i>Unused</i>	
1	B1	<i>Unused</i>	
1	B2	Scriptaid	10572
1	B4	Apicidin	10575
1	B5	HC Toxin	10576
1	B6	UNC0321 (trifluoroacetate salt)	10582
1	B7	(-)-Neplanocin A	10584
1	B8	Cl-Amidine (trifluoroacetate salt)	10599
1	B9	F-Amidine (trifluoroacetate salt)	10610
1	B10	JGB1741	10641
1	B11	coumarin-SAHA	10671
1	B12	<i>Unused</i>	
1	C1	<i>Unused</i>	
1	C2	I-BET762	10676
1	C3	UNC0638	10734
1	C4	Phthalazinone pyrazole	10735
1	C5	Isoliquiritigenin	10739
1	C6	CCG-100602	10787
1	C7	CAY10669	10974
1	C8	Zebularine	10975
1	C9	Delphinidin (chloride)	11012
1	C10	ITF 2357	11045
1	C11	UNC0631	11084
1	C12	<i>Unused</i>	
1	D1	<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, 04/10/2018

#### 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 5

Plate	Well	Contents	Item Number
1	D2	UNC0646	11085
1	D3	Methylstat (hydrate)	11091
1	D4	3-Deazaneplanocin A (hydrochloride)	11102
1	D5	Suramin (sodium salt)	11126
1	D6	Nicotinamide	11127
1	D7	2,4-Pyridinedicarboxylic Acid	11138
1	D8	PFI-1	11155
1	D9	5-Azacytidine	11164
1	D10	SGI-1027	11165
1	D11	Decitabine	11166
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	I-BET151	11181
1	E3	(+)-JQ1	11187
1	E4	(-)-JQ1	11232
1	E5	BSI-201	11304
1	E6	1-Naphthoic Acid	11322
1	E7	Sodium 4-Phenylbutyrate	11323
1	E8	Rucaparib (phosphate)	11570
1	E9	IOX1	11572
1	E10	MI-2 (hydrochloride)	11620
1	E11	MI-nc (hydrochloride)	11621
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	Gemcitabine	11690
1	F3	Lomeguatrib	11732
1	F4	GSK4112	11931
1	F5	Octyl- $\alpha$ -ketoglutarate	11970
1	F6	Daminozide	12033
1	F7	GSK-J1 (sodium salt)	12054
1	F8	GSK-J2 (sodium salt)	12056
1	F9	GSK-J4 (hydrochloride)	12073
1	F10	CI-994	12084
1	F11	CPTH2 (hydrochloride)	12086
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 5

Plate	Well	Contents	Item Number
1	G1	<i>Unused</i>	
1	G2	Etoposide	12092
1	G3	Lestaurtinib	12094
1	G4	Butyrolactone 3	12095
1	G5	Valproic Acid (sodium salt)	13033
1	G6	Tenovin-1	13085
1	G7	Tenovin-6	13086
1	G8	Sodium Butyrate	13121
1	G9	BIX01294 (hydrochloride hydrate)	13124
1	G10	Anacardic Acid	13144
1	G11	AGK2	13145
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	CAY10603	13146
1	H3	Splitomicin	13168
1	H4	CBHA	13172
1	H5	M 344	13174
1	H6	Oxamflatin	13176
1	H7	Salermide	13178
1	H8	Mirin	13208
1	H9	Pimelic Diphenylamide 106	13212
1	H10	KD 5170	13214
1	H11	Panobinostat	13280
1	H12	<i>Unused</i>	
2	A1	<i>Unused</i>	
2	A2	MS-275	13284
2	A3	HNHA	13295
2	A4	RG-108	13302
2	A5	2',3',5'-triacetyl-5-Azacytidine	13373
2	A6	S-Adenosylhomocysteine	13603
2	A7	UNC0224	13631
2	A8	Chidamide	13686
2	A9	Tubacin	13691
2	A10	3-Deazaneplanocin A	13828
2	A11	Sinefungin	13829
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 5

Plate	Well	Contents	Item Number
2	B1	<i>Unused</i>	
2	B2	Pyroxamide	13870
2	B3	N-Oxalylglycine	13944
2	B4	WDR5-0103	13945
2	B5	EPZ005687	13966
2	B6	SGC0946	13967
2	B7	UNC1215	13968
2	B8	AK-7	14004
2	B9	GSK343	14094
2	B10	Bromosporine	14119
2	B11	GSK2801	14120
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	SIRT1/2 Inhibitor IV	14407
2	C3	I-CBP112 (hydrochloride)	14468
2	C4	SGC-CBP30	14469
2	C5	UNC0642	14604
2	C6	UNC1999	14621
2	C7	(R)-PFI-2 (hydrochloride)	14678
2	C8	HPOB	15066
2	C9	2-hexyl-4-Pentynoic Acid	15205
2	C10	PFI-3	15267
2	C11	JIB-04	15338
2	C12	<i>Unused</i>	
2	D1	<i>Unused</i>	
2	D2	CAY10683	15403
2	D3	GSK126	15415
2	D4	CPI-203	15479
2	D5	6-Thioguanine	15774
2	D6	Tubastatin A	15785
2	D7	3,3'-Diindolylmethane	15927
2	D8	OTX015	15947
2	D9	5-Methylcytidine	16111
2	D10	AGK7	16152
2	D11	5-Methyl-2'-deoxycytidine	16166
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 5

Plate	Well	Contents	Item Number
2	E1	<i>Unused</i>	
2	E2	EPZ5676	16175
2	E3	MC 1568	16265
2	E4	$\alpha$ -Hydroxyglutaric Acid (sodium salt)	16374
2	E5	S-(5'-Adenosyl)-L-methionine (tosylate)	16376
2	E6	RVX-208	16424
2	E7	CUDC-101	16426
2	E8	LAQ824	16427
2	E9	Nullscript	16433
2	E10	GSK-LSD1 (hydrochloride)	16439
2	E11	RGFP966	16917
2	E12	<i>Unused</i>	
2	F1	<i>Unused</i>	
2	F2	BRD73954	16919
2	F3	<i>trans</i> -Resveratrol	70675
2	F4	DMOG	71210
2	F5	Trichostatin A	89730
2	F6	CAY10398	89740
2	F7	RSC-133	9001839
2	F8	BML-210	10005019
2	F9	Piceatannol	10009366
2	F10	CAY10591	10009797
2	F11	EX-527	10009798
2	F12	<i>Unused</i>	
2	G1	<i>Unused</i>	
2	G2	SAHA	10009929
2	G3	2-PCPA (hydrochloride)	10010494
2	G4	<i>Unused</i>	
2	G5	<i>Unused</i>	
2	G6	<i>Unused</i>	
2	G7	<i>Unused</i>	
2	G8	<i>Unused</i>	
2	G9	<i>Unused</i>	
2	G10	<i>Unused</i>	
2	G11	<i>Unused</i>	
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