

# PRODUCT INFORMATION



## Stem Cell Small Molecule Screening Library (96-Well)

Item No. 9001827 • Batch No. 0465456

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

Page 1 of 6

Plate	Well	Contents	Item Number
1	A1	<i>Unused</i>	
1	A2	Zebularine	10975
1	A3	Nicotinamide	11127
1	A4	5-Azacytidine	11164
1	A5	Sodium 4-Phenylbutyrate	11323
1	A6	Ponatinib	11494
1	A7	AT-9283	11496
1	A8	CP 690,550	11598
1	A9	Enzastaurin	11601
1	A10	Ruxolitinib	11609
1	A11	SB 505124	11793
1	A12	<i>Unused</i>	
1	B1	<i>Unused</i>	
1	B2	Kartogenin	11826
1	B3	Cyclosporin A	12088
1	B4	Cardiogenol C (hydrochloride)	13187
1	B5	Neurodazine	13224
1	B6	Gö 6983	13311
1	B7	Pifithrin- $\alpha$	13326
1	B8	Bisindolylmaleimide IX (mesylate)	13334
1	B9	LY364947	13341
1	B10	IWR-1-endo	13659
1	B11	GANT 61	13841
1	B12	<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, 04/07/2016

#### 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	C1	<i>Unused</i>	
1	C2	WIKI4	14057
1	C3	Indirubin	14155
1	C4	GANT 58	14193
1	C5	KY 02111	14315
1	C6	Thioridazine (hydrochloride)	14400
1	C7	8-bromo-Cyclic AMP (sodium salt)	14431
1	C8	1-EBIO	14575
1	C9	DBZ	14627
1	C10	TG101209	14696
1	C11	GSK3 $\beta$ Inhibitor VIII	14704
1	C12	<i>Unused</i>	
1	D1	<i>Unused</i>	
1	D2	WP1066	14736
1	D3	SANT 1	14933
1	D4	PF-05274857 (hydrochloride)	15022
1	D5	Amiodarone (hydrochloride)	15213
1	D6	ID-8	15222
1	D7	IWP-L6	15243
1	D8	AM580	15261
1	D9	GSK429286A	15262
1	D10	(+)-BAY-K-8644	15263
1	D11	LY2157299	15312
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	LY2109761	15409
1	E3	CHIR98014	15578
1	E4	PS-48	15642
1	E5	IQ-1	16153
1	E6	CH 223191	16154
1	E7	PluriSIn 1	16161
1	E8	LY411575	16162
1	E9	ISX-9	16165
1	E10	WHI-P154	16178
1	E11	BMS 833923	16240
1	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 3 of 6

Plate	Well	Contents	Item Number
1	F1	<i>Unused</i>	
1	F2	CYT387	16246
1	F3	GW 788388	16255
1	F4	ICG-001	16257
1	F5	LDE225 (phosphate)	16263
1	F6	RKI-1447	16278
1	F7	SB-525334	16281
1	F8	TG101348	16289
1	F9	XL019	16298
1	F10	Wnt-C59	16644
1	F11	AZD 1080	16676
1	F12	<i>Unused</i>	
1	G1	<i>Unused</i>	
1	G2	Astragaloside A	16677
1	G3	K02288	16678
1	G4	DMH1	16679
1	G5	P7C3	16682
1	G6	LY2784544	16705
1	G7	MK-0752	16706
1	G8	Baricitinib	16707
1	G9	SR 3677	16708
1	G10	Pacritinib	16709
1	G11	Avagacestat	16711
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	LY2811376	16712
1	H3	Semagacestat	16713
1	H4	NVP-BSK805 (hydrochloride)	16714
1	H5	Sotrastaurin	16726
1	H6	Tideglusib	16727
1	H7	AZD 2858	16728
1	H8	ZM 39923 (hydrochloride)	16729
1	H9	AZ 960	16731
1	H10	CEP-33779	16732
1	H11	1-Azakenpauillone	16733
1	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 4 of 6

Plate	Well	Contents	Item Number
2	A1	<i>Unused</i>	
2	A2	Pifithrin- $\mu$	10748
2	A3	Dexamethasone	11015
2	A4	Retinoic Acid	11017
2	A5	Forskolin	11018
2	A6	Decitabine	11166
2	A7	Cyclopamine	11321
2	A8	LDN193189	11802
2	A9	A-769662	11900
2	A10	Dorsomorphin	11967
2	A11	SB-431542	13031
2	A12	<i>Unused</i>	
2	B1	<i>Unused</i>	
2	B2	PD 173074*	13032
2	B3	Valproic Acid (sodium salt)	13033
2	B4	PD 0325901*	13034
2	B5	SB 203580	13067
2	B6	Sodium Butyrate	13121
2	B7	CHIR99021	13122
2	B8	BIO	13123
2	B9	BIX01294 (hydrochloride hydrate)	13124
2	B10	DAPT	13197
2	B11	RG-108	13302
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	IBMX	13347
2	C3	XAV939	13596
2	C4	Ku-0063794	13597
2	C5	GDC 0449	13613
2	C6	IDE1	13816
2	C7	IDE2	13817
2	C8	IWP-2	13951
2	C9	IWP-2-V2	13952
2	C10	IWP-3	13953
2	C11	IWP-4	13954
2	C12	<i>Unused</i>	

\*Sold for research purposes under agreement from Pfizer Inc.

## 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	D1	<i>Unused</i>	
2	D2	Prostaglandin E <sub>2</sub>	14010
2	D3	Thiazovivin	14245
2	D4	β-Glycerophosphate (sodium salt hydrate)	14405
2	D5	Doxycycline (hyclate)	14422
2	D6	(-)-Indolactam V	14647
2	D7	L-Ascorbic Acid	14656
2	D8	Cyclic Pifithrin-α (hydrobromide)	14748
2	D9	IPA-3	14759
2	D10	ALK5 Inhibitor II	14794
2	D11	(+)-BAY-K-8644	15263
2	D12	<i>Unused</i>	
2	E1	<i>Unused</i>	
2	E2	BI-D1870	15264
2	E3	JK 184	15265
2	E4	LY294002	70920
2	E5	Trichostatin A	89730
2	E6	Reversine	10004412
2	E7	Y-27632 (hydrochloride)	10005583
2	E8	PD 98059	10006726
2	E9	SC-1	10009557
2	E10	Purmorphamine	10009634
2	E11	Kenpaullone	10010239
2	E12	<i>Unused</i>	
2	F1	<i>Unused</i>	
2	F2	2-PCPA (hydrochloride)	10010494
2	F3	5-Methyl-2'-deoxycytidine	16166
2	F4	CD437	71610
2	F5	Troglitazone	71750
2	F6	EC 23	9002073
2	F7	Quercetin	10005169
2	F8	SAHA	10009929
2	F9	AICAR	10010241
2	F10	SB-216763	10010246
2	F11	SB-415286	10010247
2	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 6 of 6

Plate	Well	Contents	Item Number
2	G1	<i>Unused</i>	
2	G2	AG-490	10010311
2	G3	SP 600125	10010466
2	G4	HA-1077 (hydrochloride)	10010559
2	G5	TWS119	10011251
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>	
2	H1	<i>Unused</i>	
2	H2	<i>Unused</i>	
2	H3	<i>Unused</i>	
2	H4	<i>Unused</i>	
2	H5	<i>Unused</i>	
2	H6	<i>Unused</i>	
2	H7	<i>Unused</i>	
2	H8	<i>Unused</i>	
2	H9	<i>Unused</i>	
2	H10	<i>Unused</i>	
2	H11	<i>Unused</i>	
2	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