

# PRODUCT INFORMATION



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

Item No. 9001827 • Batch No. 0622233

Page 1 of 6

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

Plate	Well	Contents	Item Number	Primary Target*
1	A1	Unused		
1	A2	24(S),25-epoxy Cholesterol	10131	LXR $\alpha$
1	A3	Garcinol	10566	p300, PCAF
1	A4	StemRegenin 1	10625	AhR
1	A5	STF-31	11173	Glut1
1	A6	Ruxolitinib	11609	JAK1, JAK2
1	A7	(+)-ar-Turmerone	11759	Stem Cell Proliferation
1	A8	Kartogenin	11826	Stem Cell Differentiation
1	A9	SAG	11914	Smo
1	A10	SB-431542 (hydrate)	13031	ALK5
1	A11	PD 173074	13032	FGFR1
1	A12	Unused		
1	B1	Unused		
1	B2	PD 0325901	13034	MEK
1	B3	CHIR99021	13122	GSK3 $\alpha$ , GSK3 $\beta$
1	B4	BIO	13123	GSK3 $\alpha$ , GSK3 $\beta$
1	B5	BIX01294 (hydrochloride hydrate)	13124	G9a HMTase
1	B6	Cardiogenol C (hydrochloride)	13187	Stem Cell Differentiation
1	B7	NSC 23766 (hydrochloride)	13196	Rac1
1	B8	DAPT	13197	$\gamma$ -Secretase
1	B9	Neurodazine	13224	Stem Cell Differentiation
1	B10	VPC 23019	13240	S1P Receptors
1	B11	D 4476	13305	CK1
1	B12	Unused		
1	C1	Unused		
1	C2	EHNA (hydrochloride)	13352	Adenosine Deaminase
1	C3	Salinomycin (sodium salt)	13579	Cancer Stem Cells (CSCs)
1	C4	IDE1	13816	Stem Cell Differentiation
1	C5	IDE2	13817	Stem Cell Differentiation
1	C6	GANT 61	13841	GLI1
1	C7	IWP-2	13951	Wnt
1	C8	IWP-3	13953	Porcupine
1	C9	IWP-4	13954	Wnt
1	C10	WIKI4	14057	TNKS1, TNKS2
1	C11	GSK343	14094	EZH2
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, 08/23/2021

### 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	Primary Target*
1	D1	<i>Unused</i>		
1	D2	OAC1	14102	Oct4
1	D3	OAC2	14103	Oct4
1	D4	OAC3	14104	Oct4
1	D5	Thiazovivin	14245	ROCK
1	D6	KY 02111	14315	Stem Cell Differentiation
1	D7	Pyronin Y	14488	dsRNA
1	D8	Homoharringtonine	14631	Cancer Stem Cells (CSCs)
1	D9	16,16-dimethyl Prostaglandin E <sub>2</sub>	14750	EP Receptors
1	D10	ALK5 Inhibitor II	14794	ALK5
1	D11	ML115	15178	STAT3
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	ML-239	15188	Cancer Stem Cells (CSCs)
1	E3	ML-243	15189	Cancer Stem Cells (CSCs)
1	E4	MoTP	15204	Stem Cell Proliferaton
1	E5	Shz-1	15211	Stem Cell Differentiation
1	E6	ID-8	15222	Stem Cell Proliferaton
1	E7	IWP-L6	15243	PORCN
1	E8	AM580	15261	RAR $\alpha$
1	E9	Disulfiram	15303	Aldehyde Dehydrogenase
1	E10	LY2157299	15312	ALK1
1	E11	Galiellalactone	15387	STAT3
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	Stauprimide	15398	NME2
1	F3	GSK126	15415	EZH2
1	F4	Muscone	16067	Stem Cell Differentiation
1	F5	TCS 2210	16142	Stem Cell Differentiation
1	F6	TTNPB	16144	RARs
1	F7	IQ-1	16153	Stem Cell Pluripotency
1	F8	CH 223191	16154	AhR
1	F9	PluriSln 1	16161	Steroyl-CoA Desaturase
1	F10	LY411575	16162	$\gamma$ -Secretase
1	F11	GSA 10	16164	Smo
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	Primary Target*
1	G1	<i>Unused</i>		
1	G2	ISX-9	16165	Stem Cell Differentiation
1	G3	EPZ004777 (formate)	16173	DOT1L
1	G4	Pyrintegrin	16179	Stem Cell Survival
1	G5	NQDI-1	16226	ASK1
1	G6	BMS 833923	16240	Smo
1	G7	CYT387	16246	JAK1, JAK2
1	G8	ICG-001	16257	CBP-mediated Transcription
1	G9	LDE225 (phosphate)	16263	Smo
1	G10	PTC-209	16277	BMI-1
1	G11	TDZD-8	16287	GSK3 $\beta$
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	Dihydrolipoic Acid	16372	Stem Cell Proliferaton
1	H3	L-Ascorbic Acid 2-phosphate (magnesium salt hydrate)	16457	Stem Cell Differentiation
1	H4	Rhodamine 123 (chloride)	16672	Stem Cell Side Population
1	H5	Astragaloside A	16677	Stem Cell Differentiation
1	H6	K02288	16678	ALK1, ALK2
1	H7	DMH1	16679	ALK2
1	H8	MK-0752	16706	$\gamma$ -Secretase
1	H9	NVP-BSK805 (hydrochloride)	16714	JAK2
1	H10	Sotrastaurin	16726	PKC
1	H11	AZD 2858	16728	GSK3 $\beta$
1	H12	<i>Unused</i>		
2	A1	<i>Unused</i>		
2	A2	1-Azakenpaullone	16733	GSK3 $\beta$
2	A3	Compound 56	17254	EGFR
2	A4	Calhex 231	17268	CaSR
2	A5	Bisindolylmaleimide X (hydrochloride)	17511	PKC
2	A6	CASIN	17694	Cdc42
2	A7	SMER28	17768	Stem Cell Differentiation
2	A8	Cevimeline (hydrochloride)	17802	mAChRs
2	A9	XL147	18011	PI3K
2	A10	Ispinesib	18014	Eg5
2	A11	SW033291	18040	15-PGDH
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	Primary Target*
2	B1	<i>Unused</i>		
2	B2	A-419259 (hydrochloride)	18168	Src Family Kinases
2	B3	Hydrocortisone 21-hemisuccinate (sodium salt)	18226	Stem Cell Differentiation
2	B4	BRD-K4477	18415	Stem Cell Differentiation
2	B5	BMS 453	19076	RARs
2	B6	LLP-3	19088	Survivin
2	B7	TY 52156	19119	S1P <sub>3</sub>
2	B8	A-939572	19123	stearoyl-CoA Desaturase 1
2	B9	O4I2	19135	Oct3/4
2	B10	MS351	19549	CBX7
2	B11	Petunidin (chloride)	19755	Cancer Stem Cells (CSCs)
2	B12	<i>Unused</i>		
2	C1	<i>Unused</i>		
2	C2	Tetrandrine	19874	Stem Cells
2	C3	WZB117	19900	Glut1
2	C4	O4I1	19935	Oct3/4
2	C5	(S)-(-)-BAY-K-8644	19988	Stem Cell Reprogramming
2	C6	cis-Vaccenic Acid	20023	Stem Cell Differentiation
2	C7	BRD7116	20570	Cancer Stem Cells (CSCs)
2	C8	CP21R7	20573	GSK3 $\beta$
2	C9	AS-8351	20811	Stem Cell Reprogramming
2	C10	CGP 77675 (hydrate)	21089	Src Family Kinases
2	C11	FLI-06	21272	Notch Signaling
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	ML-390	21395	DHODH
2	D3	Bentamapimod	21482	JNKs
2	D4	Reparixin (L-Lysine salt)	21492	CXCR1, CXCR2
2	D5	BRD6125	21804	Stem Cell Differentiation
2	D6	MMP-8 Inhibitor I	21852	MMP-8
2	D7	Napabucasin	22255	Cancer Stem Cells (CSCs)
2	D8	IT-901	22298	NF- $\kappa$ B c-Rel
2	D9	Lin28 1632	22401	Lin28
2	D10	Valproic Acid Acyl-D-Glucuronide	22459	Class I HDACs
2	D11	Epiblastin A	22758	CK1 $\alpha$ , CK1 $\delta$ , CK1 $\epsilon$
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	Primary Target*
2	E1	<i>Unused</i>		
2	E2	Amcasertib	22980	Cancer Stem Cells (CSCs)
2	E3	ITD 1	23326	TGF- $\beta$
2	E4	NSC 23005 (sodium salt)	23327	p18
2	E5	FAI	23429	5S rRNA
2	E6	Hexaconazole	24118	Stem Cell Differentiation
2	E7	TA-01	24667	Stem Cell Differentiation
2	E8	TA-02	24668	Stem Cell Differentiation
2	E9	A-419259	25605	Src Family Kinases
2	E10	PD 407824	25989	Chk1, WEE1
2	E11	HhAntag	28296	Shh Signaling
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	TCS 21311	30063	JAK3
2	F3	YL-109	30290	Cancer Stem Cells (CSCs)
2	F4	PD 173955	30618	Tyrosine Kinases
2	F5	ZM 449829	32820	JAK3
2	F6	MF438	33349	stearoyl-CoA Desaturase 1
2	F7	PTC-596	33529	BMI1
2	F8	CPPG	34362	mGluRs
2	F9	GYKI 52466 (hydrochloride)	34418	AMPA Receptors
2	F10	Rosiglitazone	71740	PPAR $\gamma$
2	F11	Nanaomycin A	9001479	Stem Cell Differentiation
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	A 83-01	9001799	ALK4, ALK5, ALK7
2	G3	RSC-133	9001839	Stem Cell Reprogramming
2	G4	(R)-nitro-Blebbistatin	9001935	Unknown
2	G5	EC 23	9002073	Stem Cell Differentiation
2	G6	PF-04449913	9002936	Smo
2	G7	Reversine	10004412	Stem Cell Differentiation
2	G8	Y-27632 (hydrochloride)	10005583	ROCKs
2	G9	Retreversine	10006606	Stem Cell Differentiation
2	G10	FK-506	10007965	Calcineurin
2	G11	SC-1	10009557	Stem Cell Renewal
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 6

Plate	Well	Contents	Item Number	Primary Target*
2	H1	<i>Unused</i>		
2	H2	Purmorphamine	10009634	Stem Cell Differentiation
2	H3	Lavendustin A	10010268	EGFR-associated Tyrosine Kinase
2	H4	Tranilcypromine (hydrochloride)	10010494	LSD1
2	H5	AG-1296	10010592	PDGFR
2	H6	TWS119	10011251	GSK3 $\beta$
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>		

\*Primary Target is based on available  $IC_{50}$  and/or  $K_i$  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