

# PRODUCT INFORMATION



## Autophagy-focused Screening Library

Item No. 23537 • Batch No. 0517796

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	Unused	
1	A2	SB939	10443
1	A3	PCI 34051	10444
1	A4	Sirtinol	10523
1	A5	Scriptaid	10572
1	A6	IU1	10617
1	A7	Pifithrin- $\mu$	10748
1	A8	Piperlongumine	11006
1	A9	Delphinidin (chloride)	11012
1	A10	Forskolin	11018
1	A11	SMI-4a	11029
1	A12	Unused	
1	B1	Unused	
1	B2	ITF 2357	11045
1	B3	PFI-1	11155
1	B4	Sodium 4-Phenylbutyrate	11323
1	B5	YM-155	11490
1	B6	Obatoclox (mesylate)	11499
1	B7	ABT-737	11501
1	B8	ABT-888 (hydrochloride)	11505
1	B9	IOX2	11573
1	B10	Temsirolimus	11590
1	B11	Everolimus	11597
1	B12	Unused	
1	C1	Unused	
1	C2	AZD 1152-HQPA	11602
1	C3	Tamoxifen (citrate)	11629
1	C4	Gemcitabine	11690
1	C5	Vinblastine (sulfate)	11762
1	C6	Vincristine (sulfate)	11764
1	C7	Dioscin	11834
1	C8	Brefeldin A	11861
1	C9	A-769662	11900
1	C10	Dorsomorphin	11967
1	C11	Rottlerin	12006
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, 06/01/2023

#### 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	CI-994	12084
1	D3	2-Methoxyestradiol	13021
1	D4	Valproic Acid (sodium salt)	13033
1	D5	STF-62247	13084
1	D6	Tenovin-1	13085
1	D7	Metformin (hydrochloride)	13118
1	D8	M 344	13174
1	D9	3-Methyladenine	13242
1	D10	Tamoxifen	13258
1	D11	MS-275	13284
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	FK-866	13287
1	E3	Pifithrin- $\alpha$	13326
1	E4	Rapamycin	13346
1	E5	Geldanamycin	13355
1	E6	MK-0457	13600
1	E7	ZM 447439	13601
1	E8	MLN8237	13602
1	E9	WYE-354	13604
1	E10	PP242	13643
1	E11	Nocodazole	13857
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	JNJ-26481585 (hydrochloride)	14088
1	F3	Temozolomide	14163
1	F4	Torin 2	14185
1	F5	Chloroquine (phosphate)	14194
1	F6	2-deoxy-D-Glucose	14325
1	F7	Acridine Orange	14338
1	F8	Nimodipine	14573
1	F9	GSK2578215A	14603
1	F10	( $\pm$ )-Thalidomide	14610
1	F11	Cyclic Pifithrin- $\alpha$ (hydrobromide)	14748
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	Nafamostat (mesylate)	14837
1	G3	Amlodipine	14838
1	G4	Loperamide (hydrochloride)	14875
1	G5	Omeprazole	14880
1	G6	SNS-314 (mesylate)	14974
1	G7	Azithromycin	15004
1	G8	Doxorubicin (hydrochloride)	15007
1	G9	PF-04691502	15017
1	G10	Trifluoperazine (hydrochloride)	15068
1	G11	Amiodarone (hydrochloride)	15213
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	P005091	15224
1	H3	WP1130	15227
1	H4	Roxadustat	15294
1	H5	DBeQ	15318
1	H6	Taurolidine	15374
1	H7	LDN-57444	15396
1	H8	Carbamazepine	15422
1	H9	Ranolazine (hydrochloride)	15604
1	H10	Tubastatin A	15785
1	H11	Clonidine (hydrochloride)	15949
1	H12	<i>Unused</i>	
2	A1	<i>Unused</i>	
2	A2	<i>p</i> -nitro-Pifithrin- $\alpha$	16209
2	A3	GNE-0877	16251
2	A4	GNE-7915	16252
2	A5	JNJ-26854165	16259
2	A6	MC 1568	16265
2	A7	ONX 0912	16269
2	A8	ONX 0914	16271
2	A9	XL147 analog 2	16299
2	A10	SGI-1776	16423
2	A11	CUDC-101	16426
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	LAQ824	16427
2	B3	Bupivacaine	16618
2	B4	AZD 8055	16978
2	B5	SAR405	16979
2	B6	Perhexiline (maleate)	16982
2	B7	PIK-III	17002
2	B8	VLX600	17192
2	B9	DL- <i>threo</i> -PPMP (hydrochloride)	17236
2	B10	Verteporfin	17334
2	B11	ML-240	17373
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	OSI-027	17379
2	C3	Vps34-IN1	17392
2	C4	(±)-Felodipine	17535
2	C5	Isradipine	17536
2	C6	Nitrendipine	17549
2	C7	Carfilzomib	17554
2	C8	NMS-873	17674
2	C9	Quinacrine analog 34	17733
2	C10	CE3F4	17767
2	C11	SMER28	17768
2	C12	<i>Unused</i>	
2	D1	<i>Unused</i>	
2	D2	Spautin-1	17769
2	D3	Hydroxychloroquine (sulfate)	17911
2	D4	MHY1485	18008
2	D5	LRRK2-IN-1	18094
2	D6	CCT128930	18194
2	D7	Mocetinostat	18287
2	D8	PF-06447475	18375
2	D9	MLN9708	18386
2	D10	Danusertib	18387
2	D11	SBI-0206965	18477
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	GSK621	19155
2	E3	MK-5108	19167
2	E4	CPTH6 (hydrobromide)	19828
2	E5	Chrysophanol	19870
2	E6	MRT68921	19905
2	E7	PCI 24781	20059
2	E8	Sulfacetamide (sodium salt)	20377
2	E9	Flunarizine (hydrochloride)	20385
2	E10	L-Tetrahydropalmatine	20535
2	E11	PP121	21196
2	E12	<i>Unused</i>	
2	F1	<i>Unused</i>	
2	F2	CYC-116	21205
2	F3	Dorsomorphin (hydrochloride)	21207
2	F4	Oleuropein	21220
2	F5	Nilvadipine	21243
2	F6	Azelnidipine	21249
2	F7	ACY-1215	21531
2	F8	Aurora A Inhibitor I	21600
2	F9	ZLN024	21644
2	F10	Ro 3280	21669
2	F11	NVP-BGT226	22142
2	F12	<i>Unused</i>	
2	G1	<i>Unused</i>	
2	G2	Divalproex (sodium salt)	22960
2	G3	Clevidipine Butyrate	23025
2	G4	GNE-9605	23446
2	G5	Parthenolide	70080
2	G6	Aspirin	70260
2	G7	<i>trans</i> -Resveratrol	70675
2	G8	LY294002	70920
2	G9	Trichostatin A	89730
2	G10	(E/Z)-Belinostat	10004279
2	G11	Y-27632 (hydrochloride)	10005583
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
2	H1	<i>Unused</i>	
2	H2	RITA	10006426
2	H3	PI-103	10009209
2	H4	EX-527	10009798
2	H5	SAHA	10009929
2	H6	BAY 11-7082	10010266
2	H7	SP600125	10010466
2	H8	HA-1077 (hydrochloride)	10010559
2	H9	Wortmannin	10010591
2	H10	SRT 1720 (hydrochloride)	10011020
2	H11	(S)-MG132	10012628
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