

# PRODUCT INFORMATION



## Cell Death Screening Library

Item No. 35093 • Batch No. 0628999

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

Page 1 of 6

Plate	Well	Contents	Item Number	Major Pathway	Primary Target***
1	A1	Unused			
1	A2	Chelerythrine (chloride)*	11314	Apoptosis	PKC, Bcl-xL
1	A3	Actinomycin D*	11421	Apoptosis	DNA
1	A4	Obatoclox (mesylate)*	11499	Apoptosis	Bcl-2, Bcl-W, Bcl-xL, Mcl-1
1	A5	ABT-737*	11501	Apoptosis	Bcl-2, Bcl-W, Bcl-xL
1	A6	Camptothecin*	11694	Apoptosis	Topoisomerase I
1	A7	Cycloheximide*	14126	Apoptosis	Translation Elongation
1	A8	BAM7*	14566	Apoptosis	Bax
1	A9	(-)-Nutlin-3*	18585	Apoptosis	p53-Mdm2
1	A10	TW-37*	20999	Apoptosis	Bcl-2, Bcl-xL, Mcl-1
1	A11	A-1331852*	22963	Apoptosis	Bcl-xL
1	A12	Unused			
1	B1	Unused			
1	B2	Apoptosis Activator 2*	10004176	Apoptosis	Apaf-1
1	B3	Torin 1*	10997	Autophagy	mTORC1, mTORC2
1	B4	Dorsomorphin*	11967	Autophagy	AMPK
1	B5	Rapamycin*	13346	Autophagy	mTORC1
1	B6	Ku-0063794*	13597	Autophagy	mTORC1, mTORC2
1	B7	EN6*	28715	Autophagy	mTORC1
1	B8	BIBF 1120*	11022	Entosis	FGFR, PDGFR, VEGFR
1	B9	Artemisinin*	11816	Ferroptosis	ROS Production
1	B10	Artesunate*	11817	Ferroptosis	ROS Production
1	B11	L-Buthionine-(S,R)-Sulfoximine*	14484	Ferroptosis	γ-Glutamylcysteine Synthetase
1	B12	Unused			
1	C1	Unused			
1	C2	Auranofin*	15316	Ferroptosis	Thioredoxin Reductase
1	C3	N-Acetyl-4-benzoquinone imine*	16115	Ferroptosis	Glutathione
1	C4	Pyrimethamine*	16472	Ferroptosis	DHFR
1	C5	Erastin*	17754	Ferroptosis	System x <sub>c</sub> <sup>-</sup> Cystine-glutamate Transporter
1	C6	(1S,3R)-RSL3*	19288	Ferroptosis	GPX4
1	C7	ML-162*	20455	Ferroptosis	GPX4
1	C8	ML-390*	21395	Ferroptosis	DHODH
1	C9	ML-210*	23282	Ferroptosis	GPX4
1	C10	Pralatrexate*	23689	Ferroptosis	DHFR
1	C11	Brequinar*	24445	Ferroptosis	DHODH
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, 05/02/2023

#### 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	Major Pathway	Primary Target***
1	D1	<i>Unused</i>			
1	D2	FINO <sub>2</sub> *	25096	Ferroptosis	GPX4, Iron
1	D3	FIN56*	25180	Ferroptosis	GPX4
1	D4	Erastin2*	27087	Ferroptosis	System x <sub>c</sub> <sup>-</sup> Cystine-glutamate Transporter
1	D5	Chlorido[N,N'-disalicylidene-1,2-phenylenediamine]iron(III)*	28788	Ferroptosis	N/A
1	D6	Diaveridine*	29427	Ferroptosis	DHFR
1	D7	iFSP1*	29483	Ferroptosis	AIFM2/FSP1
1	D8	JKE-1674*	30784	Ferroptosis	GPX4
1	D9	JKE-1716*	30786	Ferroptosis	GPX4
1	D10	BAY-2402234*	33259	Ferroptosis	DHODH
1	D11	ASLAN003*	33516	Ferroptosis	DHODH
1	D12	<i>Unused</i>			
1	E1	<i>Unused</i>			
1	E2	Sorafenib*	10009644	Ferroptosis	pan-Kinase, System x <sub>c</sub> <sup>-</sup> Cystine-glutamate Transporter
1	E3	Shikonin*	14751	Necroptosis	PTEN, Pyruvate Kinase M2
1	E4	HS-173*	19156	Necroptosis	PI3Kα
1	E5	Staurosporine*	81590	Necroptosis	PKCα
1	E6	Resibufogenin*	9001348	Necroptosis	Na <sup>+</sup> /K <sup>+</sup> -ATPase
1	E7	Berberine (chloride)*	10006427	Necroptosis	N/A
1	E8	A23187*	11016	NETosis	Calcium
1	E9	Ionomycin (calcium salt)*	11932	NETosis	Calcium
1	E10	Phorbol 12-myristate 13-acetate*	10008014	NETosis	PKC
1	E11	Nigericin (sodium salt)*	11437	Pyroptosis	Potassium
1	E12	<i>Unused</i>			
1	F1	<i>Unused</i>			
1	F2	JNJ-26481585 (hydrochloride)*	14088	Pyroptosis	pan-HDAC
1	F3	<i>Unused</i>			
1	F4	<i>Unused</i>			
1	F5	<i>Unused</i>			
1	F6	<i>Unused</i>			
1	F7	Ac-DEVD-CHO (trifluoroacetate salt)**	10017	Apoptosis	pan-Caspase
1	F8	Caspase-3 Inhibitor VII**	13306	Apoptosis	Caspase-3
1	F9	PD 150606**	13859	Apoptosis	μ-Calpain, m-Calpain
1	F10	Calpain Inhibitor III**	14283	Apoptosis	μ-Calpain, m-Calpain
1	F11	Z-DEVD-FMK**	14414	Apoptosis	Caspase-3
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	Major Pathway	Primary Target***
1	G1	<i>Unused</i>			
1	G2	Z-VAD(OMe)-FMK**	14463	Apoptosis	pan-Caspase
1	G3	Caspase-3/7 Inhibitor I**	14464	Apoptosis	Caspase-3, Caspase-7
1	G4	Calpain Inhibitor I**	14921	Apoptosis	μ-Calpain, m-Calpain, Cathepsin B, Cathepsin L
1	G5	Q-VD-OPH**	15260	Apoptosis	pan-Caspase
1	G6	BI-6C9**	17265	Apoptosis	Truncated Bid
1	G7	NS 3694**	19190	Apoptosis	Apoptosome
1	G8	Emricasan**	22204	Apoptosis	pan-Caspase
1	G9	PD 151746**	23417	Apoptosis	μ-Calpain
1	G10	Caspase-9 Inhibitor III (trifluoroacetate salt)**	24306	Apoptosis	Caspase-9
1	G11	Ac-IETD-CHO (trifluoroacetate salt)**	27100	Apoptosis	Caspase-8
1	G12	<i>Unused</i>			
1	H1	<i>Unused</i>			
1	H2	CAY10746**	29814	Apoptosis	ROCK1, ROCK2
1	H3	Y-27632 (hydrochloride)**	10005583	Apoptosis	ROCK1, ROCK2, PRK2
1	H4	PD 169316**	10006727	Apoptosis	p38 MAPK
1	H5	(S)-H-1152 (hydrochloride)**	10007653	Apoptosis	ROCK
1	H6	Bafilomycin A <sub>1</sub> **	11038	Autophagy-dependent cell death	V-ATPase
1	H7	Concanamycin A**	11050	Autophagy-dependent cell death	V-ATPase
1	H8	CAY10657**	11140	Autophagy-dependent cell death	N/A
1	H9	3-Methyladenine**	13242	Autophagy-dependent cell death	PI3K
1	H10	Bafilomycin B <sub>1</sub> **	14005	Autophagy-dependent cell death	V-ATPase
1	H11	Chloroquine (phosphate)**	14194	Autophagy-dependent cell death	Autophagosome
1	H12	<i>Unused</i>			
2	A1	<i>Unused</i>			
2	A2	IMD 0354**	17290	Autophagy-dependent cell death	IKKβ
2	A3	Spautin-1**	17769	Autophagy-dependent cell death	USP10, USP13
2	A4	SBI-0206965**	18477	Autophagy-dependent cell death	ULK1, ULK2
2	A5	MRT68921**	19905	Autophagy-dependent cell death	ULK1, ULK2
2	A6	LY294002**	70920	Autophagy-dependent cell death	PI3K
2	A7	SC-514**	10010267	Autophagy-dependent cell death	IKKβ
2	A8	Wortmannin**	10010591	Autophagy-dependent cell death	PI3K, Plk1, Plk3
2	A9	ML-7 (hydrochloride)**	11801	Entotic cell death	MLCK
2	A10	(-)-Blebbistatin**	13013	Entotic cell death	Myosin IIA ATPase, Myosin IIB ATPase
2	A11	MLS-573151**	15269	Entotic cell death	Cdc42
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	Major Pathway	Primary Target***
2	B1	<i>Unused</i>			
2	B2	(S)-CCG-1423**	20107	Entotic cell death	MRTF-A
2	B3	Scriptaid**	10572	Ferroptosis	HDAC
2	B4	Rosiglitazone (maleate)**	11884	Ferroptosis	PPAR $\gamma$
2	B5	Idebenone**	15475	Ferroptosis	Free Radicals
2	B6	Ciclopirox**	16021	Ferroptosis	HIF-PH2, HIF-1 $\alpha$
2	B7	Deferasirox**	16753	Ferroptosis	Iron
2	B8	Zaragozic Acid A**	17452	Ferroptosis	Squalene Synthase
2	B9	Ferrostatin-1**	17729	Ferroptosis	ROS Production, Lipid Peroxidation
2	B10	Lipoxstatin-1**	17730	Ferroptosis	Lipid Peroxidation
2	B11	GKT137831**	17764	Ferroptosis	NOX1, NOX4
2	B12	<i>Unused</i>			
2	C1	<i>Unused</i>			
2	C2	YM-53601**	18113	Ferroptosis	Squalene Synthase
2	C3	FLT3 Inhibitor**	18462	Ferroptosis	FLT3
2	C4	Deferiprone**	20387	Ferroptosis	Iron
2	C5	SRS11-92**	25689	Ferroptosis	Iron
2	C6	Lipoxygenin**	26310	Ferroptosis	5-LO
2	C7	SRS16-86**	26752	Ferroptosis	N/A
2	C8	BLX3887**	27391	Ferroptosis	15-LO-1
2	C9	( $\pm$ )- $\alpha$ -Tocopherol Acetate**	28399	Ferroptosis	Oxidative Stress
2	C10	IM-93**	28794	Ferroptosis	Free Radicals, Lipid Peroxidation
2	C11	UAMC-3203 (hydrochloride)**	29346	Ferroptosis	N/A
2	C12	<i>Unused</i>			
2	D1	<i>Unused</i>			
2	D2	CN128**	30833	Ferroptosis	Iron
2	D3	RC363**	31407	Ferroptosis	ROS Production
2	D4	PRGL493**	32748	Ferroptosis	ACSL4
2	D5	GSK2795039**	33777	Ferroptosis	NOX2
2	D6	LY83583**	70230	Ferroptosis	Guanylate Cyclase
2	D7	Baicalein**	70610	Ferroptosis	12-LO, 15-LO-1, Lipid Peroxidation
2	D8	Troglitazone**	71750	Ferroptosis	PPAR $\gamma$
2	D9	Mitoquinol**	89950	Ferroptosis	Antioxidant
2	D10	Zileuton**	10006967	Ferroptosis	5-LO
2	D11	PD 146176**	10010518	Ferroptosis	15-LO-1
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	Major Pathway	Primary Target***
2	E1	<i>Unused</i>			
2	E2	Deferoxamine (mesylate)**	14595	Lysosome-dependent cell death	Prolyl Hydroxylase
2	E3	Cathepsin G Inhibitor I**	14928	Lysosome-dependent cell death	Cathepsin G
2	E4	CA-074 methyl ester**	18469	Lysosome-dependent cell death	Cathepsin B, Cathepsin L
2	E5	Odanacatib**	21466	Lysosome-dependent cell death	Cathepsin K
2	E6	Cathepsin L Inhibitor**	23249	Lysosome-dependent cell death	Cathepsin L
2	E7	Necrostatin-5**	10527	Necroptosis	RIPK1
2	E8	Necrostatin-7**	10528	Necroptosis	N/A
2	E9	Necrostatin-2**	11657	Necroptosis	N/A
2	E10	Necrostatin-1**	11658	Necroptosis	RIPK1
2	E11	Dabrafenib**	16989	Necroptosis	pan-Kinase
2	E12	<i>Unused</i>			
2	F1	<i>Unused</i>			
2	F2	Necrox-5 (methanesulfonate)**	17278	Necroptosis	ROS Production, RNS Production
2	F3	Necrosulfonamide**	20844	Necroptosis	MLKL
2	F4	GSK872**	23300	Necroptosis	RIPK3
2	F5	HS-1371**	31033	Necroptosis	RIPK3
2	F6	Cl-Amidine (hydrochloride)**	10599	NETotic cell death	PAD1, PAD3, PAD4
2	F7	4-Aminobenzoic Acid hydrazide**	14845	NETotic cell death	MPO
2	F8	GSK484 (hydrochloride)**	17488	NETotic cell death	PAD4
2	F9	Sivelestat (sodium salt hydrate)**	17779	NETotic cell death	Neutrophil Elastase
2	F10	GW 311616A**	27957	NETotic cell death	Neutrophil Elastase
2	F11	AZD 7986**	33184	NETotic cell death	DPP-1
2	F12	<i>Unused</i>			
2	G1	<i>Unused</i>			
2	G2	N-acetyl-L-Carnosine**	18817	Oxeiptosis	Lipid Peroxidation
2	G3	BSI-201**	11304	Parthanatos	PARP1
2	G4	1,5-Isoquinolinediol**	14438	Parthanatos	PARP1
2	G5	NMS-P118**	21340	Parthanatos	PARP1
2	G6	BYK204165**	21430	Parthanatos	PARP1
2	G7	AG-14361**	24677	Parthanatos	PARP1
2	G8	Ac-YVAD-CMK**	10014	Pyroptosis	Caspase-1
2	G9	Ac-YVAD-CHO**	10016	Pyroptosis	Caspase-1
2	G10	Wedelolactone**	11796	Pyroptosis	NF-κB
2	G11	Glyburide (potassium salt)**	15009	Pyroptosis	SUR1/K <sub>ir</sub> 6.2
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	Major Pathway	Primary Target***
2	H1	<i>Unused</i>			
2	H2	Disulfiram**	15303	Pyroptosis	Aldehyde Dehydrogenase
2	H3	MCC950**	17510	Pyroptosis	NLRP3
2	H4	Dapansutrine**	24671	Pyroptosis	NLRP3
2	H5	Ac-LEVD-CHO (trifluoroacetate salt)**	27433	Pyroptosis	Caspase-4
2	H6	INF39**	28476	Pyroptosis	NLRP3
2	H7	CY-09**	33353	Pyroptosis	NLRP3
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>			

\* Inducer of the cell death pathway

\*\* Inhibitor of the cell death pathway

\*\*\* 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