

PRODUCT INFORMATION



Immuno-Oncology Screening Library

Item No. 28760 • Batch No. 0581972

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

Page 1 of 4

Plate	Well	Contents	Item Number
1	A1	<i>Unused</i>	
1	A2	PLX4032	10618
1	A3	I-BET762	10676
1	A4	IC-87114	11589
1	A5	AMD 3465 (hydrobromide)	11959
1	A6	INCB 3284 (dimesylate)	11963
1	A7	SB 297006	11964
1	A8	SR 3335	12072
1	A9	Pazopanib	12097
1	A10	CAY10603	13146
1	A11	Sunitinib (malate)	13159
1	A12	<i>Unused</i>	
1	B1	<i>Unused</i>	
1	B2	MS-275	13284
1	B3	SB 225002	13336
1	B4	WZ811	13639
1	B5	HDAC-IN-7	13686
1	B6	Axitinib	13813
1	B7	TAK-242	13871
1	B8	Chloroquine (phosphate)	14194
1	B9	DMXAA	14617
1	B10	R-848	14806
1	B11	Imiquimod	14956
1	B12	<i>Unused</i>	
1	C1	<i>Unused</i>	
1	C2	CAL-101	15279
1	C3	LY2157299	15312
1	C4	OTX015	15947
1	C5	Ibrutinib	16274
1	C6	Trametinib	16292
1	C7	SR 1078	16503
1	C8	SD 208	16619
1	C9	Dabrafenib	16989
1	C10	Encorafenib	16994
1	C11	Binimetinib	16996
1	C12	<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, 04/21/2021

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 4

Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	CGS 21680 (hydrochloride)	17126
1	D3	RS 504393	17330
1	D4	GSK2126458	17377
1	D5	GSK2269557	17388
1	D6	MRS1706	17417
1	D7	Defactinib	17737
1	D8	Hydroxychloroquine (sulfate)	17911
1	D9	Dyphylline	17958
1	D10	AVL-292	17993
1	D11	RS 102895 (hydrochloride)	18207
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	INCB024360 analog	18265
1	E3	Mocetinostat	18287
1	E4	CGI1746	18414
1	E5	BX 471	18503
1	E6	Doxofylline	18746
1	E7	Regadenoson	18814
1	E8	Gardiquimod	18877
1	E9	PD-1/PD-L1 Inhibitor 2	19132
1	E10	Lenvatinib	19375
1	E11	IDO-IN-1	19402
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	SR 0987	19503
1	F3	Cobimetinib	19563
1	F4	GS-9620	19628
1	F5	Epacadostat	19875
1	F6	Acalabrutinib	19899
1	F7	PD-1/PD-L1 Inhibitor 1	19914
1	F8	Vipadenant	20239
1	F9	ZM 241385	20447
1	F10	Ticlopidine (hydrochloride)	20770
1	F11	Proxiphylline	20937
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 4

Plate	Well	Contents	Item Number
1	G1	<i>Unused</i>	
1	G2	GSK2981278	20974
1	G3	Danirixin	20980
1	G4	5'-N-Ethylcarboxamidoadenosine (hydrate)	21420
1	G5	Reparixin (L-Lysine salt)	21492
1	G6	(±)-NLG-919	21509
1	G7	Theobromine	21745
1	G8	MK-0812	21803
1	G9	Sulfamethoxypyridazine	21875
1	G10	BEC (hydrochloride)	22145
1	G11	Preladenant	22210
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	G10	22353
1	H3	Istradefylline	22958
1	H4	Vercirnon	23131
1	H5	Sulfamethizole	23355
1	H6	MSX-122	23426
1	H7	Theophylline	23760
1	H8	EOS200271	25003
1	H9	CPI-444	25024
1	H10	Atractylenolide I	25061
1	H11	Baohuoside I	25062
1	H12	<i>Unused</i>	
2	A1	<i>Unused</i>	
2	A2	CCCP	25458
2	A3	H-151	25857
2	A4	C-176	25859
2	A5	C-178	25860
2	A6	AZD 4635	26322
2	A7	IPI-549	26416
2	A8	TAK-779	26744
2	A9	C29	27029
2	A10	2-Cl-IB-MECA	27336
2	A11	STING Agonist C11	27886
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 4

Plate	Well	Contents	Item Number
2	B1	<i>Unused</i>	
2	B2	AZD 5069	28297
2	B3	GSK805	9002444
2	B4	ABH	10006862
2	B5	Sorafenib	10009644
2	B6	SAHA	10009929
2	B7	Plerixafor (hydrochloride hydrate)	10011332
2	B8	<i>Unused</i>	
2	B9	<i>Unused</i>	
2	B10	<i>Unused</i>	
2	B11	<i>Unused</i>	
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	<i>Unused</i>	
2	C3	<i>Unused</i>	
2	C4	<i>Unused</i>	
2	C5	<i>Unused</i>	
2	C6	<i>Unused</i>	
2	C7	<i>Unused</i>	
2	C8	<i>Unused</i>	
2	C9	<i>Unused</i>	
2	C10	<i>Unused</i>	
2	C11	<i>Unused</i>	
2	C12	<i>Unused</i>	
2	D1	<i>Unused</i>	
2	D2	<i>Unused</i>	
2	D3	<i>Unused</i>	
2	D4	<i>Unused</i>	
2	D5	<i>Unused</i>	
2	D6	<i>Unused</i>	
2	D7	<i>Unused</i>	
2	D8	<i>Unused</i>	
2	D9	<i>Unused</i>	
2	D10	<i>Unused</i>	
2	D11	<i>Unused</i>	
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