

PRODUCT INFORMATION



Comprehensive Kinase Screening Library

Item No. 28940 • Batch No. 0582322

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

Page 1 of 30

Plate	Well	Contents	Item Number	Primary Target*
1	A1	Unused		
1	A2	Necrostatin-5	10527	RIPK2
1	A3	MK-8669	10543	mTOR
1	A4	AZD 1480	10702	JAK2
1	A5	SMI-4a	11029	PIMs
1	A6	BAY 61-3606 (hydrochloride)	11423	Syk
1	A7	CHR-6494	11478	Haspin
1	A8	Cediranib	11495	VEGFR1, VEGFR2, VEGFR3
1	A9	Saracatinib	11497	c-Src, Abl
1	A10	GSK1059615	11569	PI3K α
1	A11	CI-1040	11580	MEK
1	A12	Unused		
1	B1	Unused		
1	B2	IC-87114	11589	PI3K p110 δ
1	B3	MK-2206 (hydrochloride)	11593	Akt1, Akt2, Akt3
1	B4	AZD 6244	11599	MEK1, MEK2
1	B5	Enzastaurin	11601	PKC family
1	B6	K145 (hydrochloride)	11691	SPHK2
1	B7	Embelin	11838	XIAP
1	B8	AZD 4547	13111	FGFR1, FGFR2, FGFR3
1	B9	Gö 6976	13310	PKC α , PKC β 1
1	B10	IKK-16 (hydrochloride)	13313	IKK α , IKK β , IKK complex
1	B11	PD 158780	13329	HER2, HER3, HER4
1	B12	Unused		
1	C1	Unused		
1	C2	1-NM-PP1	13330	Analog-sensitive mutant kinases
1	C3	SB 218078	13335	Chk1
1	C4	ST638	13337	Tyrosine kinases
1	C5	CAY10621	13371	SPHK1
1	C6	Arcyriaflavin A	13516	Cdk4, CaMKII
1	C7	VEGFR2 Kinase Inhibitor I	13578	VEGFR2
1	C8	WYE-354	13604	mTOR
1	C9	NVP-231	13858	Ceramide kinase
1	C10	LY333531 (hydrochloride)	13964	PKC β 1, PKC β 2
1	C11	AD57 (hydrochloride)	13975	RET
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, 09/04/2024

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