

PRODUCT INFORMATION



Kinase Screening Library (96-Well)

Item No. 10505 • Batch No. 0668513

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

Page 1 of 6

Plate	Well	Contents	Item Number	Primary Target*
1	A1	Unused		
1	A2	TG003	10398	CLK1, CLK4
1	A3	Doramapimod	10460	p38 MAPK
1	A4	Erlotinib	10483	EGFR
1	A5	PLX4032	10618	B-RAF, B-RAV600E
1	A6	Phthalazinone pyrazole	10735	Aurora A
1	A7	AG-879	10793	BMX
1	A8	Torin 1	10997	mTOR
1	A9	BIBF 1120	11022	FGFR, PDGFR, VEGFR
1	A10	SMI-4a	11029	PIMs
1	A11	AS-703026	11226	MEK1, MEK2
1	A12	Unused		
1	B1	Unused		
1	B2	Chelerythrine (chloride)	11314	Bcl-xL, PKC
1	B3	R406	11422	Syk
1	B4	AZD 7762	11491	Chk1, Chk2
1	B5	Dasatinib	11498	Src, Abl, Lck
1	B6	GSK1059615	11569	PI3K α
1	B7	Ruxolitinib	11609	JAK1, JAK2
1	B8	Necrostatin-1	11658	RIPK1
1	B9	INK128	11811	mTOR
1	B10	Bosutinib	12030	c-Src, Abl
1	B11	Canertinib (hydrochloride)	12076	EGFR
1	B12	Unused		
1	C1	Unused		
1	C2	(R)-Crizotinib	12087	c-Met, ALK
1	C3	SB-431542 (hydrate)	13031	ALK5
1	C4	PD 173074*	13032	FGFR1
1	C5	PD 0325901*	13034	MEK
1	C6	SB 203580	13067	p38 MAPK
1	C7	VX-702	13108	p38 MAPK
1	C8	Emodin	13109	CK2
1	C9	CHIR99021	13122	GSK3
1	C10	BIO	13123	GSK3
1	C11	Imatinib (mesylate)	13139	Multi-kinase
1	C12	Unused		

*Sold for research purposes under agreement from Pfizer Inc.

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, 03/01/2024

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

PRODUCT INFORMATION



Page 2 of 6

Plate	Well	Contents	Item Number	Primary Target*
1	D1	<i>Unused</i>		
1	D2	Sunitinib (malate)	13159	FLK1, FLT3, PDGFR β
1	D3	PP2	13198	Fyn, HCK, LYN
1	D4	3-Methyladenine	13242	PI3K
1	D5	Bisindolylmaleimide I	13298	PKC
1	D6	Bisindolylmaleimide V	13300	S6K
1	D7	D 4476	13305	CK1
1	D8	NU 7026	13308	DNA-PK
1	D9	Gö 6983	13311	PKC
1	D10	H-9 (hydrochloride)	13312	PKG
1	D11	Indirubin-3'-monoxime	13314	GSK3 β
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	NU 6102	13317	Cdk1, Cdk2
1	E3	KN-62	13318	CAMKII
1	E4	KN-93	13319	CAMKII
1	E5	CGP 57380	13322	MNK1
1	E6	(S)-Glycyl-H-1152 (hydrochloride)	13332	ROCK-II
1	E7	Bisindolylmaleimide IX (mesylate)	13334	GSK3, PKC
1	E8	SU6656	13338	Yes
1	E9	LY364947	13341	ALK5
1	E10	CAY10621	13371	SPHK1
1	E11	YM-201636	13576	PIKfyve
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	ZM 447439	13601	Aurora B
1	F3	AS-041164	13622	PI3K γ
1	F4	NVP-AEW541 (hydrochloride)	13641	IGF-1R
1	F5	PP242	13643	mTOR
1	F6	ABT-869	13653	VEGFR family, PDGFR family
1	F7	CAY10622	13687	ROCK-I, ROCK-II
1	F8	CAY10626	13838	PI3K α , mTOR
1	F9	SU 6668	13873	Aurora B
1	F10	BMS-536924	14666	IGF-1R, InsR
1	F11	PHA-665752	14703	c-Met
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	BX-912	14708	PDK1
1	G3	Vatalanib (hydrochloride)	14868	VEGFR family
1	G4	KRN 633	14871	VEGFR family
1	G5	LCK Inhibitor	15135	LCK
1	G6	BI-D1870	15264	RSKs
1	G7	CRT0066101 (hydrochloride)	15337	PKD
1	G8	GNE-7915	16252	LRRK2
1	G9	GNF-5	16254	Bcr-Abl
1	G10	BMS 345541 (hydrochloride)	16667	IKK α , IKK β
1	G11	NVP-TAE226	17685	FAK, PYK2 β
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	AZ191	17693	DYRK1B
1	H3	Tie2 Kinase Inhibitor	17858	Tie2
1	H4	Afuresertib (hydrochloride)	17988	PKB/Akt
1	H5	DDR1-IN-1 (hydrate)	18092	DDR1
1	H6	GSK2334470	18095	PDK1
1	H7	JNK Inhibitor XVI	18096	JNKs
1	H8	BI-6727	18193	Plks
1	H9	Syk Inhibitor II	18202	Syk
1	H10	PHA-767491 (hydrochloride)	18218	Cdc7
1	H11	PF-06463922	18371	ALK
1	H12	<i>Unused</i>		
2	A1	<i>Unused</i>		
2	A2	PF-562271 (besylate)	18499	FAK
2	A3	BMS-777607	18517	Axl family, Met family
2	A4	URMC-099	19147	LRRK2, MLKs
2	A5	BGJ398	19157	FGFR family
2	A6	Mps1-IN-1	19175	Mps1
2	A7	WZ4003	19177	NUAK1, NUAK2
2	A8	MLCK Inhibitor Peptide 18	19181	MLCK
2	A9	BIX 02565	19183	RSK2
2	A10	AP26113	19778	ALK
2	A11	AZ 3146	19991	Mps1
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	AZD 1208	20235	PIMs
2	B3	RPI-1	20384	RET
2	B4	LDN-211904	20453	EphB3
2	B5	BMS-5	21072	LIMK1, LIMK2
2	B6	TAS 120	21136	FGFR family
2	B7	Apatinib	21268	VEGFR2
2	B8	ARQ 092	21388	PKB/Akt
2	B9	LY2606368	21490	Chk1
2	B10	FRAX597	22205	PAK1, PAK2, PAK3
2	B11	KW 2449	22207	FLT3, Abl
2	B12	<i>Unused</i>		
2	C1	<i>Unused</i>		
2	C2	UNC569	22966	Axl family
2	C3	XMD16-5	23353	TNK2
2	C4	Tilfrinib	25647	BRK/PTK6
2	C5	AZD 0156	25648	ATM
2	C6	LOXO-101	27056	Trk family
2	C7	LOXO-195	27062	Trk family
2	C8	U-0126	70970	MEK, AMPK
2	C9	Staurosporine	81590	PKC
2	C10	Y-27632 (hydrochloride)	10005583	PRK2, ROCK1, ROCK2
2	C11	Leelamine	10006148	PDK
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	PD 169316	10006727	p38 MAPK
2	D3	TGX-221	10007349	PI3K p110 β
2	D4	(S)-H-1152 (hydrochloride)	10007653	ROCK
2	D5	AS-605240	10007707	PI3K γ
2	D6	JNJ-10198409	10008131	PDGFR
2	D7	CAY10505	10009078	CK2
2	D8	PI-103	10009209	DNA-PK, PI3K, mTOR
2	D9	PIK-75 (hydrochloride)	10009210	PI3K p110 α
2	D10	Sphingosine Kinase Inhibitor 2	10009222	SPHK1
2	D11	SC-1	10009557	RasGAP, ERK1
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	(R)-Roscovitine	10009569	CDKs
2	E3	Sorafenib	10009644	pan-Kinase, System xc-Cystine-glutamate Transporter
2	E4	CAY10561	10010043	ERK2
2	E5	PI3-Kinase α Inhibitor 2	10010177	PI3K α
2	E6	ML-9	10010236	Multi-kinase
2	E7	Erbstatin analog	10010238	EGFR
2	E8	Kenpaullone	10010239	CDKs, GSK3 β
2	E9	Olomoucine	10010240	CDKs
2	E10	AG-494	10010242	EGFR
2	E11	AG-825	10010243	ErbB2
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	AG-1478	10010244	EGFR
2	F3	SB-216763	10010246	GSK3
2	F4	H-8 (hydrochloride)	10010249	PKA, PKG
2	F5	LFM-A13	10010265	BTK
2	F6	SC-514	10010267	IKK β
2	F7	RG-13022	10010309	EGFR
2	F8	AG-490	10010311	JAK2
2	F9	AG-183	10010315	EGFR
2	F10	Lavendustin C	10010329	EGFR
2	F11	GW 5074	10010368	Raf-1
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	5-Iodotubercidin	10010375	CK1, ERK2, PKC
2	G3	SB 202190	10010399	p38 MAPK
2	G4	CAY10571	10010400	p38 α MAPK
2	G5	Nilotinib	10010422	Bcr-Abl
2	G6	SP600125	10010466	JNKs
2	G7	H-89 (hydrochloride)	10010556	Multi-kinase
2	G8	HA-1077 (hydrochloride)	10010559	ROCK-II
2	G9	AG-370	10010568	PDGFR
2	G10	Wortmannin	10010591	PI3K, PIk1, PIk3
2	G11	AG-1296	10010592	PDGFR
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	WHI-P131	10011246	JAK3
2	H3	CAY10574	10011247	Cdk9
2	H4	CAY10576	10011249	IKK ϵ
2	H5	NH125	10011250	EEF2K
2	H6	TWS119	10011251	GSK3 β
2	H7	CAY10578	10011264	CK2
2	H8	PD 184161	10012431	MEK1, MEK2
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