

PRODUCT INFORMATION



Kinase Screening Library (96-Well)

Item No. 10505 • Batch No. 0612092

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	Doramapimod	10460	p38 MAPK
1	A3	Erlotinib	10483	EGFR
1	A4	Torin 1	10997	mTOR
1	A5	AZD 7762	11491	Chk1, Chk2
1	A6	Dasatinib	11498	Src
1	A7	GSK1059615	11569	PI3K α
1	A8	Ruxolitinib	11609	JAK1, JAK2
1	A9	Necrostatin-1	11658	RIP1
1	A10	INK128	11811	mTOR
1	A11	Bosutinib	12030	c-Src, Abl
1	A12	Unused		
1	B1	Unused		
1	B2	(R)-Crizotinib	12087	c-Met, ALK5
1	B3	SB-431542 (hydrate)	13031	ALK5
1	B4	PD 173074*	13032	FGFR1
1	B5	PD 0325901*	13034	MEK
1	B6	SB 203580	13067	p38 MAPK
1	B7	Emodin	13109	CK2
1	B8	CHIR99021	13122	GSK3
1	B9	BIO	13123	GSK3
1	B10	Imatinib (mesylate)	13139	PDGFR, c-Kit
1	B11	3-Methyladenine	13242	PI3K
1	B12	Unused		
1	C1	Unused		
1	C2	Bisindolylmaleimide V	13300	S6K
1	C3	D 4476	13305	CK1
1	C4	NU 6102	13317	Cdk1, Cdk2
1	C5	Gö 6983	13311	PKC
1	C6	Indirubin-3'-monoxime	13314	GSK3 β
1	C7	NU 6102	13308	DNA-PK
1	C8	KN-62	13318	CAMKII
1	C9	SU6656	13338	Yes
1	C10	YM-201636	13576	PIKfyve

*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, 05/13/2021

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 6

Plate	Well	Contents	Item Number	Target
1	C11	ZM 447439	13601	Aurora B
1	C12	<i>Unused</i>		
1	D1	<i>Unused</i>		
1	D2	PP242	13643	mTOR
1	D3	ABT-869	13653	VEGFR family, PDGFR family
1	D4	CAY10626	13838	PI3K α , mTOR
1	D5	PHA-665752	14703	c-Met
1	D6	LCK Inhibitor	15135	LcK
1	D7	BI-D1870	15264	RSKs
1	D8	GNE-7915	16252	LRRK2
1	D9	BMS 345541 (trifluoroacetate salt)	16667	IKK α , IKK β
1	D10	NVP-TAE226	17685	FAK, PYK2 β
1	D11	Tie2 Kinase Inhibitor	17858	Tie2
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	BI-6727	18193	Plks
1	E3	Syk Inhibitor II	18202	Syk
1	E4	PHA-767491 (hydrochloride)	18218	Cdc7
1	E5	PF-06463922	18371	ALK
1	E6	PF-562271 (besylate)	18499	FAK
1	E7	Mps1-IN-1	19175	Mps1
1	E8	MLCK Inhibitor Peptide 18	19181	MLCK
1	E9	TAS 120	21136	FGFR family
1	E10	Apatinib	21268	VEGFR2
1	E11	ARQ-092	21388	PKB/Akt
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	LY2606368	21490	Chk1
1	F3	FRAX597	22205	PAK1, PAK2, PAK3
1	F4	KW 2449	22207	FLT3, Abl
1	F5	<i>Unused</i>		
1	F6	LOXO-195	27062	Trk family
1	F7	U-0126	70970	MEK1, MEK2
1	F8	Staurosporine	81590	PKC
1	F9	PD 169316	10006727	p38 MAPK
1	F10	AS-605240	10007707	PI3K γ

*Sold for research purposes under agreement from Pfizer Inc.

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	Target
1	F11	PI-103	10009209	DNA-PK, PI3K, mTOR
1	F12	<i>Unused</i>		
1	G1	<i>Unused</i>		
1	G2	Sphingosine Kinase Inhibitor 2	10009222	SPHK1
1	G3	SC-1	10009557	ERK1
1	G4	(R)-Roscovitine	10009569	CDKs
1	G5	ML-9	10010236	Multi-kinase
1	G6	Olomoucine	10010240	CDKs
1	G7	AG-494	10010242	EGFR
1	G8	AG-825	10010243	ErbB2
1	G9	SB-216763	10010246	GSK3
1	G10	AG-490	10010311	JAK2
1	G11	AG-183	10010315	EGFR
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	Lavendustin C	10010329	EGFR
1	H3	SB 202190	10010399	p38 MAPK
1	H4	CAY10571	10010400	p38 α MAPK
1	H5	Nilotinib	10010422	Bcr-Abl
1	H6	SP 600125	10010466	JNK1, JNK2, JNK3, JNKs
1	H7	H-89 (hydrochloride)	10010556	Multi-kinase
1	H8	AG-1296	10010592	PDGFR
1	H9	WHI-P131	10011246	JAK3
1	H10	CAY10576	10011249	IKK ϵ
1	H11	TWS119	10011251	GSK3 β
1	H12	<i>Unused</i>		
2	A1	<i>Unused</i>		
2	A2	TG003	10398	CLK1, CLK4
2	A3	PLX4032	10618	B-RAF
2	A4	Phthalazinone pyrazole	10735	Aurora A
2	A5	AG-879	10793	BMX
2	A6	BIBF 1120	11022	Multi-kinase
2	A7	SMI-4a	11029	PIMs
2	A8	AS-703026	11226	MEK1, MEK2
2	A9	Chelerythrine (chloride)	11314	PKC
2	A10	R406	11422	Syk

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	Target
2	A11	Canertinib (hydrochloride)	12076	EGFR
2	A12	<i>Unused</i>		
2	B1	<i>Unused</i>		
2	B2	VX-702	13108	p38 MAPK
2	B3	Sunitinib (malate)	13159	FLK1, PDGFR β
2	B4	Bisindolylmaleimide I	13298	PKC
2	B5	H-9 (hydrochloride)	13312	PKG
2	B6	KN-93	13319	CAMKII
2	B7	CGP 57380	13322	MNK1
2	B8	(S)-Glycyl-H-1152 (hydrochloride)	13332	ROCK-II
2	B9	Bisindolylmaleimide IX (mesylate)	13334	GSK3, PKC
2	B10	LY364947	13341	ALK5
2	B11	CAY10621	13371	SPHK1
2	B12	<i>Unused</i>		
2	C1	<i>Unused</i>		
2	C2	AS-041164	13622	PI3Ky
2	C3	NVP-AEW541 (hydrochloride)	13641	IGF-1R
2	C4	CAY10622	13687	ROCK-I, ROCK-II
2	C5	17 β -hydroxy Wortmannin	13812	PI3K
2	C6	SU 6668	13873	Aurora B
2	C7	BMS-536924	14666	IGF-1R, InsR
2	C8	BX-912	14708	PDK1
2	C9	Vatalanib (hydrochloride)	14868	VEGFR family
2	C10	KRN 633	14871	VEGFR family
2	C11	CRT0066101 (hydrochloride)	15337	PKD
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	GNF-5	16254	Bcr-Abl
2	D3	AZ191	17693	DYRK1B
2	D4	Afuresertib (hydrochloride)	17988	PKB/Akt
2	D5	GSK2334470	18095	PDK1
2	D6	JNK Inhibitor XVI	18096	JNK1, JNK2, JNK3, JNKs
2	D7	BMS-777607	18517	Axl family, Met family
2	D8	URMC-099	19147	LRRK2, MLKs
2	D9	BGJ398	19157	FGFR family
2	D10	WZ4003	19177	NUAK1, NUAKE2

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	Target
2	D11	AP26113	19778	ALK
2	D12	<i>Unused</i>		
2	E1	<i>Unused</i>		
2	E2	AZ 3146	19991	Mps1
2	E3	AZD 1208	20235	PIMs
2	E4	RPI-1	20384	RET
2	E5	LDN-211904	20453	EphB3
2	E6	BMS-5	21072	LIMK1, LIMK2
2	E7	UNC569	22966	Axl family
2	E8	XMD16-5	23353	TNK2
2	E9	Tilfrinib	25647	BRK/PTK6
2	E10	AZD 0156	25648	ATM
2	E11	LOXO-101	27056	Trk family
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	Y-27632 (hydrochloride)	10005583	ROCK-I
2	F3	Leelamine	10006148	PDK
2	F4	TGX-221	10007349	PI3K p110 β
2	F5	(S)-H-1152 (hydrochloride)	10007653	ROCK
2	F6	JNJ-10198409	10008131	PDGFR
2	F7	CAY10505	10009078	CK2
2	F8	PIK-75 (hydrochloride)	10009210	PI3K p110 α
2	F9	Sorafenib	10009644	Raf-1
2	F10	CAY10561	10010043	ERK2
2	F11	PI3-Kinase α Inhibitor 2	10010177	PI3K α
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	Erbstatin analog	10010238	EGFR
2	G3	Kenpaullone	10010239	CDKs, GSK3 β
2	G4	AG-1478	10010244	EGFR
2	G5	H-8 (hydrochloride)	10010249	PKA, PKG
2	G6	LFM-A13	10010265	BTK
2	G7	SC-514	10010267	IKK2
2	G8	RG-13022	10010309	EGFR
2	G9	GW 5074	10010368	Raf-1

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	Target
2	G10	5-Iodotubercidin	10010375	CK1, ERK2, PKC
2	G11	HA-1077 (hydrochloride)	10010559	ROCK-II
2	G12	<i>Unused</i>		
2	H1	<i>Unused</i>		
2	H2	AG-370	10010568	PDGFR
2	H3	Wortmannin	10010591	PI3K
2	H4	CAY10574	10011247	Cdk9
2	H5	NH125	10011250	EEF2K
2	H6	CAY10578	10011264	CK2
2	H7	PD 184161	10012431	MEK1, MEK2
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>		

*Sold for research purposes under agreement from Pfizer Inc.

**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