

PRODUCT INFORMATION



Kinase Screening Library (96-Well)

Item No. 10505 • Batch No. 0521277

The composition of the library will always vary somewhat depending upon our inventory.

Page 1 of 6

Plate	Well	Contents	Item Number	Activity
1	A1	Unused		
1	A2	TG003	10398	Clk1/4 Inhibitor
1	A3	PKC 412	10459	PKC inhibitor
1	A4	Doramapimod	10460	MAPK inhibitor
1	A5	Paclitaxel	10461	Kinase inhibitor
1	A6	Erlotinib	10483	EGFR inhibitor
1	A7	NVP-BEZ235	10565	PI3K/mTOR inhibitor
1	A8	Phthalazinone pyrazole	10735	Aurora-A kinase
1	A9	AG-879	10793	Tyrosine kinase inhibitor
1	A10	1-NA-PP1	10954	Src family kinase inhibitor
1	A11	Torin 1	10997	mTOR inhibitor
1	A12	Unused		
1	B1	Unused		
1	B2	BIBF 1120	11022	VEGFR inhibitor
1	B3	SMI-4a	11029	Pim-kinase inhibitor
1	B4	CAY10657	11140	IKK2 inhibitor
1	B5	AS-703026	11226	MEK inhibitor
1	B6	Chelerythrine (chloride)	11314	PKC inhibitor
1	B7	Tunicamycin	11445	EGFR/ERB inhibitor
1	B8	AZD 7762	11491	Chk1/2 inhibitor
1	B9	GSK1059615	11569	PI3K α inhibitor
1	B10	Ruxolitinib	11609	JAK1/2 Inhibitor
1	B11	Necrostatin-1	11658	RIP1 inhibitor
1	B12	Unused		
1	C1	Unused		
1	C2	SB-505124	11793	ALK inhibitor
1	C3	INK128	11811	TORC1/2 inhibitor
1	C4	Canertinib (hydrochloride)	12076	Irreversible HER family RTK inhibitor
1	C5	SB-431542	13031	TGF β /ALK inhibitor
1	C6	PD 173074*	13032	FGFR inhibitor
1	C7	Valproic Acid (sodium salt)	13033	GSK3 β inhibitor
1	C8	PD 0325901*	13034	MEK inhibitor
1	C9	SB 203580	13067	p38 MAPK inhibitor
1	C10	VX-702	13108	MAPK inhibitor

*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, 04/24/2018

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	Activity
1	C11	Emodin	13109	Casein kinase II inhibitor
1	C12	<i>Unused</i>		
1	D1	<i>Unused</i>		
1	D2	CHIR99021	13122	GSK3 α / β inhibitor
1	D3	BIO	13123	GSK3 α / β inhibitor
1	D4	Imatinib (mesylate)	13139	Tyrosine kinase inhibitor
1	D5	Sunitinib (malate)*	13159	Tyrosine kinase inhibitor
1	D6	Gefitinib	13166	EGFR inhibitor
1	D7	PP2	13198	Src family kinase inhibitor
1	D8	3-Methyladenine	13242	PI3K inhibitor
1	D9	Bisindolylmaleimide I	13298	PKC inhibitor
1	D10	Bisindolylmaleimide IV	13299	PKC inhibitor
1	D11	Bisindolylmaleimide V	13300	S6K inhibitor
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	NSC 663284	13303	Cdc25 Phosphatase Inhibitor II
1	E3	D 4476	13305	Casein kinase inhibitor
1	E4	NU 7026	13308	DNA dependent protein kinase inhibitor
1	E5	Gö 6983	13311	PKC inhibitor
1	E6	H-9 (hydrochloride)	13312	PK inhibitor
1	E7	Indirubin-3'-monoxime	13314	GSK3 β inhibitor
1	E8	NU 6102	13317	Cdk inhibitor
1	E9	KN-62	13318	CaMKII inhibitor
1	E10	KN-93	13319	CaMKII inhibitor
1	E11	CGP 57380	13322	MNK1 inhibitor
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	<i>Iso</i> -Olomoucine	13325	Inactive isomer of olomoucine
1	F3	(S)-Glycyl-H-1152 (hydrochloride)	13332	ROCK inhibitor
1	F4	Bisindolylmaleimide VIII (acetate)	13333	PKC inhibitor
1	F5	Bisindolylmaleimide IX (mesylate)	13334	PKC inhibitor
1	F6	ST638	13337	Tyrosine kinase inhibitor
1	F7	SU6656	13338	Src family kinase inhibitor
1	F8	LY364947	13341	TGF- β RI inhibitor
1	F9	SB 203580 (hydrochloride)	13344	PDK/MAPK inhibitor
1	F10	CAY10621	13371	SPHK 1 inhibitor 5C

*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	Activity
1	F11	YM-201636	13576	PIKfyve inhibitor
1	F12	<i>Unused</i>		
1	G1	<i>Unused</i>		
1	G2	ZM 447439	13601	Aurora kinase B inhibitor
1	G3	AS-041164	13622	PI3K inhibitor
1	G4	NVP-AEW541 (hydrochloride)	13641	IGF-1R Kinase inhibitor
1	G5	PP242	13643	mTORC inhibitor
1	G6	ABT-869	13653	VEGFR/RTK inhibitor
1	G7	CAY10622	13687	ROCK inhibitor
1	G8	17 β -hydroxy Wortmannin	13812	PI3K inhibitor
1	G9	CAY10626	13838	PI3K α /mTOR inhibitor
1	G10	SU 6668	13873	RTK inhibitor
1	G11	PHA-767491	18218	Cdc7 kinase inhibitor
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	N,N-Dimethylsphingosine	62575	SPHK inhibitor
1	H3	LY294002	70920	PI3K inhibitor
1	H4	U-0126	70970	MEK1/2 inhibitor
1	H5	Staurosporine	81590	PKC inhibitor
1	H6	KN-92 (hydrochloride)	9000890	CaMKII inhibitor
1	H7	AS-605240 (potassium salt)	9000980	PI3K γ inhibitor
1	H8	PD 166326	9000988	RTK/Src inhibitor
1	H9	O-1918	10004914	Selective antagonist of abnormal cannabidiol
1	H10	Y-27632 (hydrochloride)	10005583	ROCK inhibitor
1	H11	Leelamine	10006148	PDK inhibitor
1	H12	<i>Unused</i>		
2	A1	<i>Unused</i>		
2	A2	PD 98059	10006726	MAPK pathway inhibitor
2	A3	PD 169316	10006727	p38 MAPK inhibitor
2	A4	TGX-221	10007349	PI3K p110 β inhibitor
2	A5	(S)-H-1152 (hydrochloride)	10007653	ROCK inhibitor
2	A6	AS-605240	10007707	PI3K γ inhibitor
2	A7	D-erythro-Sphingosine C-18	10007907	PKC/PAP inhibitor
2	A8	JNJ-10198409	10008131	PDGF tyrosine kinase inhibitor
2	A9	Leelamine (hydrochloride)	10008614	PDK inhibitor
2	A10	Arachidonic Acid Leelamide	10008617	Fatty acid amide

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	Activity
2	A11	Lauric Acid Leelamide	10008618	Fatty acid amide
2	A12	<i>Unused</i>		
2	B1	<i>Unused</i>		
2	B2	AS-252424	10009052	PI3Ky inhibitor
2	B3	CAY10505	10009078	PI3K Inhibitor
2	B4	PI-103	10009209	PI3K inhibitor
2	B5	PIK-75 (hydrochloride)	10009210	PI3K p110 α inhibitor
2	B6	Sphingosine Kinase Inhibitor 2	10009222	SPHK 1 inhibitor
2	B7	Piceatannol	10009366	A resveratrol analog/serine/threonine protein kinase inhibitor
2	B8	SC-1	10009557	ERK1/Ras-GAP inhibitor
2	B9	(R)-Roscovitine	10009569	Cdk2/cyclin E inhibitor
2	B10	BAY 43-9006	10009644	B-Raf inhibitor
2	B11	CAY10561	10010043	ERK2 inhibitor
2	B12	<i>Unused</i>		
2	C1	<i>Unused</i>		
2	C2	AS-604850	10010175	PI3Ky inhibitor
2	C3	PI3-Kinase α Inhibitor 2	10010177	PI3K α inhibitor
2	C4	CAY10567	10010233	Akt translocation inhibitor
2	C5	ML-9	10010236	PKB/Akt inhibitor
2	C6	Triciribine	10010237	Akt activation inhibitor
2	C7	Erbstatin Analog	10010238	EGFR tyrosine kinase inhibitor
2	C8	Kenpauillone	10010239	Cdk inhibitor
2	C9	Olomoucine	10010240	Cdk inhibitor
2	C10	AG-494	10010242	EGF receptor kinase inhibitor
2	C11	AG-825	10010243	Her2/Neu tyrosine kinase activity inhibitor
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	AG-1478	10010244	EGFR kinase inhibitor
2	D3	SB-216763	10010246	GSK3 α inhibitor
2	D4	SB-415286	10010247	GSK3 α inhibitor
2	D5	AG-17	10010248	EGF receptor kinase inhibitor
2	D6	H-8 (hydrochloride)	10010249	PKA and PKG inhibitor
2	D7	LFM-A13	10010265	BTK and Plk inhibitor
2	D8	SC-514	10010267	IKK2 inhibitor
2	D9	Apigenin	10010275	Casein kinase II inhibitor
2	D10	AG-18	10010300	EGF receptor kinase inhibitor

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	Activity
2	D11	DRB	10010302	CTD kinase inhibitor
2	D12	<i>Unused</i>		
2	E1	<i>Unused</i>		
2	E2	RG-13022	10010309	EGF receptor kinase inhibitor
2	E3	RG-14620	10010310	EGF receptor kinase inhibitor
2	E4	AG-490	10010311	JAK2 activity inhibitor
2	E5	AG-82	10010312	EGF receptor kinase inhibitor
2	E6	AG-99	10010313	EGF receptor kinase inhibitor
2	E7	AG-213	10010314	EGF receptor kinase inhibitor
2	E8	AG-183	10010315	EGF receptor kinase inhibitor
2	E9	Lavendustin C	10010329	EGF receptor kinase inhibitor
2	E10	ZM 336372	10010367	ATP-competitive inhibitor of Raf-1
2	E11	5-Iodotubercidin	10010375	Broad kinase inhibitor
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	SB 202190	10010399	MAPK inhibitor
2	F3	CAY10571	10010400	Analog of SB 203580, inhibits IL-1 production
2	F4	Nilotinib	10010422	BCR/Abl kinase inhibitor
2	F5	SP 600125	10010466	JNK inhibitor
2	F6	L-threo-Sphingosine C-18	10010541	A D-erythro-sphingosine analog
2	F7	H-89 (hydrochloride)	10010556	PKA inhibitor
2	F8	HA-1077 (hydrochloride)	10010559	ROCK inhibitor, potent vasodilator
2	F9	AG-370	10010568	PDGF receptor kinase inhibitor
2	F10	Wortmannin	10010591	PI3K inhibitor
2	F11	AG-1296	10010592	PDGF receptor kinase inhibitor
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	Janex 1	10011246	JAK3 inhibitor
2	G3	CAY10574	10011247	Cdk9 inhibitor
2	G4	CAY10575	10011248	Benzimidazole analog, IKK-ε inhibitor
2	G5	CAY10576	10011249	Benzimidazole analog, IKK-ε inhibitor
2	G6	NH125	10011250	An antibacterial imidazole
2	G7	TWS119	10011251	GSK3β inhibitor
2	G8	NSC 210902	10011255	CK2 inhibitor
2	G9	CAY10577	10011256	CK2 inhibitor

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	Activity
2	G10	CAY10578	10011264	CK2 inhibitor
2	G11	PD 184161*	10012431	MEK1/2 inhibitor
2	G12	<i>Unused</i>		
2	H1	<i>Unused</i>		
2	H2	CCT018159	10012591	Hsp90 inhibitor
2	H3	Myricetin	10012600	A potent antioxidant
2	H4	<i>Unused</i>		
2	H5	<i>Unused</i>		
2	H6	<i>Unused</i>		
2	H7	<i>Unused</i>		
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.

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