

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, 06/01/2023

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 30

Plate	Well	Contents	Item Number	Primary Target*
1	D1	<i>Unused</i>		
1	D2	SD169	14156	p38α MAPK, p38β MAPK
1	D3	Amlexanox	14181	IKKε, TBK1
1	D4	PP1 (Src Inhibitor)	14244	Src family
1	D5	MS-1020	14273	JAK3
1	D6	3CAI	14303	Akt1, Akt2
1	D7	MK2 Inhibitor IV	14399	MK2
1	D8	Alsterpaullone	14428	CDKs, GSK3α, GSK3β
1	D9	Borrelidin	14436	CDK
1	D10	FAK Inhibitor 14	14485	FAK1
1	D11	AG-1295	14529	PDGFR
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	Src Kinase Inhibitor I	14592	Src family, RIPK2
1	E3	BX-912	14708	PDK1
1	E4	Shikonin	14751	PKM2
1	E5	W-7 (hydrochloride)	14826	CaMK
1	E6	Syk Inhibitor	14829	Syk
1	E7	TBCA	14831	CK2
1	E8	SU11274	14861	c-Met
1	E9	PS-1145	14862	IKKβ
1	E10	Akt Inhibitor VIII	14870	Akt1, Akt2
1	E11	NU 7441	14881	DNA-PK
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	GNF-5837	14883	Trk family
1	F3	Arctigenin	14913	Akt
1	F4	PF-573228	14924	FAK
1	F5	BX-795	14932	Multi-kinase
1	F6	HA-100 (hydrochloride)	14955	PKA, PKC, PKG
1	F7	TAS 301	14976	CaMKII
1	F8	TPCA-1	15115	IKKβ
1	F9	PLX4720	15142	B-RAF
1	F10	JAK Inhibitor I	15146	JAK family
1	F11	Cdk2 Inhibitor II	15154	Cdk2
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 30

Plate	Well	Contents	Item Number	Primary Target*
1	G1	<i>Unused</i>		
1	G2	MNS	15206	Syk
1	G3	AZD 6482	15250	PI3K β
1	G4	CAL-101	15279	PI3K p110 δ
1	G5	LY2584702 (tosylate)	15320	S6K1
1	G6	PKR Inhibitor	15323	PKR
1	G7	STO-609 (acetate)	15325	CaMKK α , CaMKK β
1	G8	YZ9	15352	PFKFB
1	G9	EGFR Inhibitor	15363	EGFR
1	G10	Butein	15371	Src
1	G11	GW 2580	15411	c-FMS
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	HNMPA	15543	InsR
1	H3	FR180204	15544	ERK1, ERK2
1	H4	ERK Inhibitor	15944	ERK
1	H5	TIC10	16109	Akt, ERK1, ERK2
1	H6	NQDI-1	16226	ASK1
1	H7	Foretinib	16249	PDGFR family, VEGFR family
1	H8	GNE-0877	16251	LRRK2
1	H9	SRPIN340	16284	SRPK1
1	H10	BIO-Acetoxime	16329	GSK3 α , GSK3 β
1	H11	Ku-55933	16336	ATM
1	H12	<i>Unused</i>		
2	A1	<i>Unused</i>		
2	A2	TG003	10398	CLK1, CLK4
2	A3	Erlotinib	10483	EGFR, Cdk2
2	A4	PLX4032	10618	B-RAF
2	A5	Torin 1	10997	mTOR
2	A6	AS-703026	11226	MEK1, MEK2
2	A7	Lapatinib	11493	EGFR, HER2
2	A8	Ponatinib	11494	Bcr-Abl
2	A9	AT-9283	11496	Multi-kinase
2	A10	Dasatinib	11498	Src family
2	A11	Temsirolimus	11590	mTOR
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 30

Plate	Well	Contents	Item Number	Primary Target*
2	B1	<i>Unused</i>		
2	B2	Tofacitinib (citrate)	11598	JAK1, JAK2, JAK3
2	B3	AZD 1152-HQPA	11602	Aurora B
2	B4	Ruxolitinib	11609	JAK1, JAK2
2	B5	α -acetyl Boswellic Acid	11673	IKKs
2	B6	LDN-193189	11802	ALK1, ALK2, ALK3, ALK6
2	B7	Dorsomorphin	11967	AMPK
2	B8	Pazopanib	12097	VEGFR family, PDGFR family
2	B9	VX-702	13108	p38 α MAPK, p38 β MAPK
2	B10	CHIR99021	13122	GSK3 α , GSK3 β
2	B11	3-Methyladenine	13242	PI3K
2	B12	<i>Unused</i>		
2	C1	<i>Unused</i>		
2	C2	Bisindolylmaleimide IV	13299	PKC
2	C3	Gö 6983	13311	PKC
2	C4	Indirubin-3'-monoxime	13314	GSK3 β
2	C5	KN-62	13318	CaMKII
2	C6	Bisindolylmaleimide IX (mesylate)	13334	PKC, GSK3
2	C7	LY364947	13341	ALK5
2	C8	Torin 2	14185	mTOR
2	C9	Chloroquine (phosphate)	14194	Non-kinase inhibitor control
2	C10	GSK3 β Inhibitor VIII	14704	GSK3 β
2	C11	ALK5 Inhibitor II	14794	ALK5
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	LY2109761	15409	ALKs
2	D3	CHIR98014	15578	GSK3 α , GSK3 β
2	D4	GW 788388	16255	ALK5
2	D5	PD 0332991 (hydrochloride)	16273	Cdk4, Cdk6
2	D6	Ibrutinib	16274	BTK
2	D7	BMS 345541 (trifluoroacetate salt)	16667	IKK β
2	D8	K02288	16678	ALK1, ALK2, ALK3, ALK6
2	D9	DMH1	16679	ALK2
2	D10	Pacritinib	16709	FLT3, JAK2, TYK2
2	D11	Sotrastaurin	16726	PKC
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 30

Plate	Well	Contents	Item Number	Primary Target*
2	E1	<i>Unused</i>		
2	E2	Tideglusib	16727	GSK3 β
2	E3	CEP-33779	16732	JAK2
2	E4	PIK-III	17002	VPS34
2	E5	BI-2536	17385	Plk1
2	E6	Vps34-IN1	17392	VPS34
2	E7	kb NB 142-70	18002	PKD
2	E8	LRRK2-IN-1	18094	LRRK2
2	E9	PHA-767491	18218	Cdc7, Cdk9
2	E10	GDC-0980	18303	PI3K, mTOR
2	E11	XL184	18464	Multi-kinase
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	Regorafenib	18498	Multi-kinase
2	F3	MK-5108	19167	Aurora A
2	F4	Lenvatinib	19375	VEGFR family, PDGFR family
2	F5	MRT68921	19905	ULK1, ULK2
2	F6	MK-1775	21266	WEE1
2	F7	CNX-774	21475	BTK
2	F8	<i>trans</i> -Resveratrol	70675	Multi-kinase
2	F9	U-0126	70970	MEK1, MEK2
2	F10	PD 169316	10006727	p38 MAPK
2	F11	TGX-221	10007349	PI3K p110 β
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	(S)-H-1152 (hydrochloride)	10007653	ROCK
2	G3	AS-252424	10009052	PI3K γ
2	G4	Piceatannol	10009366	Multi-kinase
2	G5	SC-1	10009557	ERK1
2	G6	(R)-Roscovitine	10009569	CDKs
2	G7	Sorafenib	10009644	Raf-1
2	G8	Kenpaullone	10010239	GSK3 β and CDKs
2	G9	Olomoucine	10010240	Cdk5/p35
2	G10	Apigenin	10010275	CK2
2	G11	AG-18	10010300	EGFR
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 30

Plate	Well	Contents	Item Number	Primary Target*
2	H1	Unused		
2	H2	AG-490	10010311	JAK2
2	H3	AG-82	10010312	EGFR
2	H4	ZM 336372	10010367	Raf-1
2	H5	SB 202190	10010399	p38 α MAPK, p38 β MAPK
2	H6	CAY10571	10010400	CSAID binding protein/MAPK14
2	H7	Nilotinib	10010422	Bcr-Abl
2	H8	SP600125	10010466	JNKs
2	H9	HA-1077 (hydrochloride)	10010559	ROCK-II
2	H10	AG-370	10010568	PDGFR
2	H11	CAY10575	10011248	IKK ϵ
2	H12	Unused		
3	A1	Unused		
3	A2	NVP-BEZ235	10565	PI3K, mTOR
3	A3	Phthalazinone pyrazole	10735	Aurora A
3	A4	AG-879	10793	BMX
3	A5	1-NA-PP1	10954	Analog-sensitive mutant kinases
3	A6	Chelerythrine (chloride)	11314	PKC
3	A7	Necrostatin-1	11658	RIPK1
3	A8	INK128	11811	mTOR
3	A9	Canertinib (hydrochloride)	12076	EGFR, HER4
3	A10	Lestaurtinib	12094	JAK2, PRK1
3	A11	SB-431542 (hydrate)	13031	ALK5, ALK4
3	A12	Unused		
3	B1	Unused		
3	B2	PD 173074	13032	FGFR1
3	B3	PD 0325901	13034	MEK
3	B4	Masitinib	13105	c-Kit
3	B5	Emodin	13109	CK2
3	B6	Imatinib (mesylate)	13139	PDGFR, c-Kit
3	B7	Sunitinib (malate)	13159	FLK1, PDGFR β
3	B8	Gefitinib	13166	EGFR
3	B9	PP2	13198	LCK, Fyn, HCK
3	B10	D 4476	13305	CK1
3	B11	NU 7026	13308	DNA-PK
3	B12	Unused		

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 7 of 30

Plate	Well	Contents	Item Number	Primary Target*
3	C1	Unused		
3	C2	H-9 (hydrochloride)	13312	PKG
3	C3	KN-93	13319	CaMKII
3	C4	CGP 57380	13322	MNK1
3	C5	Iso-Olomoucine	13325	Negative control
3	C6	(S)-Glycyl-H-1152 (hydrochloride)	13332	ROCK-II
3	C7	Bisindolylmaleimide VIII (acetate)	13333	PKC
3	C8	SU6656	13338	Src family
3	C9	Ku-0063794	13597	mTORC1, mTORC2
3	C10	MK-0457	13600	Aurora kinases
3	C11	MLN8237	13602	Aurora A
3	C12	Unused		
3	D1	Unused		
3	D2	AS-041164	13622	PI3Ky
3	D3	PP242	13643	mTOR
3	D4	ABT-869	13653	VEGFR family, PDGFR family
3	D5	CAY10622	13687	ROCK-I, ROCK-II
3	D6	Axitinib	13813	VEGFR family, PDGFR family
3	D7	CAY10626	13838	PI3K α , mTOR
3	D8	SU 6668	13873	Aurora B, Aurora C, PDGFR β
3	D9	Indirubin	14155	GSK3
3	D10	Thiazovivin	14245	ROCK
3	D11	PHA-665752	14703	c-Met
3	D12	Unused		
3	E1	Unused		
3	E2	Vatalanib (hydrochloride)	14868	VEGFR family
3	E3	GSK429286A	15262	ROCK
3	E4	GNE-7915	16252	LRRK2
3	E5	Trametinib	16292	MEK1, MEK2
3	E6	ZM 39923 (hydrochloride)	16729	JAK3
3	E7	Dabrafenib	16989	B-RAF, B-RAF ^{V600E} , C-RAF
3	E8	GSK2126458	17377	PI3K
3	E9	Afuresertib (hydrochloride)	17988	Akt1
3	E10	JNK Inhibitor XVI	18096	JNKs
3	E11	PF-06463922	18371	ALK, ROS1
3	E12	Unused		

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 8 of 30

Plate	Well	Contents	Item Number	Primary Target*
3	F1	Unused		
3	F2	SBI-0206965	18477	ULK1
3	F3	PF-562271 (besylate)	18499	FAK, PYK2
3	F4	Ceritinib	19374	ALK
3	F5	AZ 3146	19991	Mps1
3	F6	AZD 1208	20235	PIMs
3	F7	RPI-1	20384	RET
3	F8	BMS-5	21072	LIMK1, LIMK2
3	F9	NVP-BGT226	22142	PI3K
3	F10	FRAX597	22205	PAK1, PAK2, PAK3
3	F11	Unused		
3	F12	Unused		
3	G1	Unused		
3	G2	N,N-Dimethylsphingosine (d18:1)	62575	SPHK
3	G3	LY294002	70920	PI3K
3	G4	PD 166326	9000988	Bcr-Abl, c-Src
3	G5	Y-27632 (hydrochloride)	10005583	ROCK-I
3	G6	Sphingosine (d18:1)	10007907	PKC
3	G7	JNJ-10198409	10008131	PDGFR β
3	G8	PI-103	10009209	DNA-PK, PI3K, mTOR
3	G9	CAY10561	10010043	ERK2
3	G10	AS-604850	10010175	PI3K γ
3	G11	ML-9	10010236	Multi-kinase
3	G12	Unused		
3	H1	Unused		
3	H2	AG-1478	10010244	EGFR
3	H3	LFM-A13	10010265	BTK
3	H4	DRB	10010302	Cdk9, CK2
3	H5	AG-99	10010313	EGFR
3	H6	AG-183	10010315	EGFR
3	H7	CAY10574	10011247	Cdk9
3	H8	TWS119	10011251	GSK3 β
3	H9	CAY10578	10011264	CK2
3	H10	PD 184161	10012431	MEK1, MEK2
3	H11	Myricetin	10012600	MEK1
3	H12	Unused		

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 9 of 30

Plate	Well	Contents	Item Number	Primary Target*
4	A1	<i>Unused</i>		
4	A2	Doramapimod	10460	p38 MAPK
4	A3	BIBF 1120	11022	Multi-kinase
4	A4	Suramin (sodium salt)	11126	PKC
4	A5	R406	11422	Syk
4	A6	Tunicamycin Mixture	11445	EGFR family, IGF-1R
4	A7	AZD 7762	11491	Chk1, Chk2
4	A8	Everolimus	11597	mTOR
4	A9	Rottlerin	12006	PKC δ , MAPKAPK-5
4	A10	Bosutinib	12030	c-Src, Abl, EphB2
4	A11	BIO	13123	GSK3
4	A12	<i>Unused</i>		
4	B1	<i>Unused</i>		
4	B2	Bisindolylmaleimide I	13298	PKC
4	B3	NU 6102	13317	Cdk2, Cdk1
4	B4	Rapamycin	13346	mTORC1
4	B5	YM-201636	13576	PIKfyve
4	B6	ZM 447439	13601	Aurora B
4	B7	KRN 633	14871	VEGFR family
4	B8	Lck Inhibitor	15135	LCK
4	B9	R 1530	15255	Multi-kinase
4	B10	LY2157299	15312	ALK5/TGF β R1 kinase
4	B11	CRT0066101 (hydrochloride)	15337	PKD
4	B12	<i>Unused</i>		
4	C1	<i>Unused</i>		
4	C2	Ku-0060648	15344	DNA-PK, PI3K α , PI3K β , PI3K δ
4	C3	Xanthohumol	15399	PDPK1
4	C4	AZD 5363	15406	Akt kinases, S6K1, PKA
4	C5	ON-01910 (sodium salt)	15553	Plk1
4	C6	Fialuridine	15867	HSV thymidine kinases
4	C7	JNK Inhibitor VIII	15946	JNK1, JNK2
4	C8	CGK733	16242	ATR, ATM
4	C9	SB-525334	16281	ALK5/TGF β R1 kinase
4	C10	TG101348	16289	JAK2, FLT3, RET
4	C11	VS-5584	16294	mTOR, PI3K
4	C12	<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 10 of 30

Plate	Well	Contents	Item Number	Primary Target*
4	D1	<i>Unused</i>		
4	D2	CUDC-101	16426	EGFR, HER2
4	D3	A-443654	16499	Akt kinases
4	D4	D-Mannoheptulose	16548	Glucokinases, hexokinases
4	D5	UNC2025	16613	Mer, FLT3
4	D6	SD 208	16619	ALK5/TGFβR1 kinase
4	D7	SB-590885	16643	B-RAF, C-RAF
4	D8	AZD 1080	16676	GSK3
4	D9	Baricitinib	16707	JAK1, JAK2
4	D10	SR 3677	16708	ROCK-II
4	D11	AZD 2858	16728	GSK3β
4	D12	<i>Unused</i>		
4	E1	<i>Unused</i>		
4	E2	AZ 960	16731	JAK2
4	E3	IPI-145	16800	PI3Kγ
4	E4	OTSSP167 (hydrochloride)	16873	MELK
4	E5	GSK690693	16891	Multi-kinase
4	E6	JX-401	16898	p38α MAPK
4	E7	GW 441756	16915	TrkA
4	E8	GDC-0879	16993	B-RAF ^{V600E}
4	E9	Ku-60019	17502	ATM
4	E10	TAE684	17670	ALK
4	E11	NVP-TAE226	17685	FAK, PYK2-β
4	E12	<i>Unused</i>		
4	F1	<i>Unused</i>		
4	F2	GSK2334470	18095	PDK1
4	F3	SCH 900776	18131	Chk1
4	F4	BI-6727	18193	Plk1, Plk2
4	F5	Menaquinone 4	18423	IKK
4	F6	URMC-099	19147	LRRK2, MLKs
4	F7	BGJ398	19157	FGFR family
4	F8	SCH 772984	19166	ERK1, ERK2
4	F9	AMG 900	19176	Aurora kinases
4	F10	LDN-211904	20453	EphB3
4	F11	Emetine (hydrochloride hydrate)	21048	IkBa phosphorylation
4	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 11 of 30

Plate	Well	Contents	Item Number	Primary Target*
4	G1	<i>Unused</i>		
4	G2	TAS 120	21136	FGFR family
4	G3	Apatinib	21268	VEGFR2
4	G4	KW 2449	22207	FLT3, Abl, Abl ^{T315I}
4	G5	BLU-285	22985	KIT ^{D816V} , PDGFR α ^{D842V}
4	G6	XMD16-5	23353	TNK2
4	G7	Tilfrinib	25647	BRK/PTK6
4	G8	AZD 0156	25648	ATM
4	G9	LOXO-195	27062	Trk family, including TrkA and TrkC mutants
4	G10	Reversine	10004412	Mps1
4	G11	AS-605240	10007707	PI3K γ , PI3K α
4	G12	<i>Unused</i>		
4	H1	<i>Unused</i>		
4	H2	PIK-75 (hydrochloride)	10009210	PI3K p110 α
4	H3	Erbstatin analog	10010238	EGFR
4	H4	AG-494	10010242	EGFR
4	H5	SB-216763	10010246	GSK3
4	H6	SC-514	10010267	IKK β
4	H7	RG-13022	10010309	EGFR
4	H8	GW 5074	10010368	Raf-1
4	H9	H-89 (hydrochloride)	10010556	Multi-kinase
4	H10	Wortmannin	10010591	PI3K
4	H11	AG-1296	10010592	PDGFR
4	H12	<i>Unused</i>		
5	A1	<i>Unused</i>		
5	A2	TCS PIM-1 1	16286	Pim-1
5	A3	AZD 8055	16978	mTOR
5	A4	SAR405	16979	VPS34
5	A5	Rp-Cyclic AMPS (triethylammonium salt)	16985	PKA I , PKA II
5	A6	BYL719	16986	PI3K α
5	A7	Encorafenib	16994	B-RAF ^{V600E}
5	A8	Refametinib	16997	MEK1, MEK2
5	A9	TAK-733	16998	MEK1
5	A10	GSK650394	17001	SGK1
5	A11	PF-543 (hydrochloride)	17034	SPHK1
5	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 12 of 30

Plate	Well	Contents	Item Number	Primary Target*
5	B1	Unused		
5	B2	CH5183284	17120	FGFR1, FGFR2, FGFR3
5	B3	IC261	17260	CK1 δ , CK1 ϵ
5	B4	IMD 0354	17290	IKK β
5	B5	Picropodophyllotoxin	17329	IGF-1R
5	B6	Rhein	17345	Non-kinase inhibitor control
5	B7	GSK2606414	17376	PERK
5	B8	AZD 2014	17378	mTORC1, mTORC2
5	B9	GSK2636771	17380	PI3K p110 β
5	B10	ZSTK474	17381	PI3K p110 β , PI3K p110 γ , PI3K p110 δ
5	B11	AZD 8186	17384	PI3K β , PI3K δ
5	B12	Unused		
5	C1	Unused		
5	C2	GDC-0032	17387	PI3K α , PI3K δ , PI3K γ
5	C3	SAR260301	17391	PI3K p110 β
5	C4	SU 5614	17410	PDGFR family, VEGFR family
5	C5	(5Z)-7-Oxozeaenol	17459	TAK-1
5	C6	BMS 754807	17505	IGF-1R, InsR
5	C7	Bisindolylmaleimide X (hydrochloride)	17511	PKC
5	C8	Aurora Kinase Inhibitor II	17541	Aurora A
5	C9	JNK Inhibitor V	17542	JNK3
5	C10	DMPQ (hydrochloride)	17551	PDGFR β
5	C11	Chk2 Inhibitor II	17552	CK2
5	C12	Unused		
5	D1	Unused		
5	D2	EGFR/ErbB2 Inhibitor	17568	EGFR, HER2
5	D3	VE-821	17587	ATR
5	D4	PD 174265	17649	EGFR
5	D5	BIIB-057	17652	Syk
5	D6	LEE011	17666	Cdk4, Cdk6
5	D7	Filgotinib	17669	JAK1, JAK2
5	D8	PFK15	17689	PFKFB3
5	D9	Linsitinib	17708	IGF-1R, InsR
5	D10	Toceranib	17714	Multi-kinase
5	D11	L-779,450	17807	B-RAF
5	D12	Unused		

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 13 of 30

Plate	Well	Contents	Item Number	Primary Target*
5	E1	<i>Unused</i>		
5	E2	Pim-1 Inhibitor 2	17809	Pim-1
5	E3	3MB-PP1	17860	Plk1
5	E4	R547	17905	Cdk1, Cdk2, Cdk4
5	E5	6H05 (trifluoroacetate salt)	17985	K-Ras ^{G12C}
5	E6	Quizartinib	17986	FLT3
5	E7	CZC-54252	17998	LRRK2
5	E8	LDN-214117	18006	ALK1, ALK2
5	E9	PLX647	18012	FMS, Kit
5	E10	VX-745	18075	p38 α MAPK
5	E11	PD 153035	18080	EGFR
5	E12	<i>Unused</i>		
5	F1	<i>Unused</i>		
5	F2	SGX523	18093	c-Met
5	F3	A-419259 (hydrochloride)	18168	Src family
5	F4	Pexidartinib	18271	CSF1R/FMS, c-Kit
5	F5	MCB-613	18315	Non-kinase inhibitor control
5	F6	WYE-125132	18372	mTOR
5	F7	Danuseritib	18387	Multi-kinase
5	F8	PHA-793887	18388	CDKs
5	F9	Neratinib	18404	HER2
5	F10	AEE788	18416	EGFR family, VEGFR family
5	F11	CH5424802	18516	ALK
5	F12	<i>Unused</i>		
5	G1	<i>Unused</i>		
5	G2	Safingol	18624	PKC, SPHK
5	G3	RWJ 67657	18718	p38 α MAPK, p38 β MAPK
5	G4	Oclacitinib	18722	JAK family
5	G5	ALK-IN-1	18775	ALK
5	G6	CEP-32496	18776	B-RAF ^{V600E}
5	G7	Syk Inhibitor II (hydrochloride)	18805	Syk
5	G8	CC-401	18867	JNK1, JNK2, JNK3
5	G9	Crenolanib	18873	PDGFR family
5	G10	BAW 2881	18910	VEGFR2
5	G11	CL 387,785	19081	EGFR
5	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 14 of 30

Plate	Well	Contents	Item Number	Primary Target*
5	H1	Unused		
5	H2	KJ Pyr 9	19116	c-Myc
5	H3	SB 239063	19142	p38 MAPK
5	H4	KIRA6	19151	IRE1α
5	H5	HS-173	19156	PI3Kα
5	H6	TX1-85-1	19164	HER3
5	H7	BMX-IN-1	19165	BMX, BTK
5	H8	BMS-265246	19168	Cdk1, Cdk2
5	H9	Bafetinib	19169	Bcr-Abl
5	H10	SBE 13 (hydrochloride)	21241	Plk1
5	H11	AMG 337	21333	c-Met
5	H12	Unused		
6	A1	Unused		
6	A2	AAL-993	16351	VEGFR family
6	A3	BMS-3	19421	LIMK1, LIMK2
6	A4	(S)-p38 MAPK Inhibitor III	19696	p38 MAPK
6	A5	NSC 109555	19811	Chk2, BRK
6	A6	LJH685	19913	RSK1, RSK2, RSK3
6	A7	LJI308	19924	RSK1, RSK2, RSK3
6	A8	MGCD-265	20097	Multi-kinase
6	A9	SU 4312	20213	VEGFR2, PDGFR
6	A10	PH-797804	20244	p38α MAPK
6	A11	XL147 analog	20245	PI3Ks
6	A12	Unused		
6	B1	Unused		
6	B2	RN-486	20536	BTK
6	B3	PQ401	20979	IGF-1R
6	B4	CVT-313	21044	Cdk2
6	B5	Bohemine	21073	CDKs
6	B6	CEP-37440	21278	FAK1, ALK
6	B7	PIK-293	21283	PI3K p110δ
6	B8	Ensartinib	21341	ALK
6	B9	Aminopurvalanol A	21424	CDKs
6	B10	KX2-391	21429	Src
6	B11	H-7 (hydrochloride)	21434	PKC, PKA, PKG, PKM
6	B12	Unused		

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 15 of 30

Plate	Well	Contents	Item Number	Primary Target*
6	C1	<i>Unused</i>		
6	C2	Amuvatinib	21461	c-Kit, PDGFR α , c-MET
6	C3	Rebastinib	21465	Abl1, Abl1 ^{T315I}
6	C4	PHA-848125	21474	Cdk2/cyclin A
6	C5	Bentamapimod	21482	JNKs
6	C6	Gilteritinib	21503	FLT3, Axl
6	C7	LY3009120	21504	A-RAF, B-RAF, C-RAF
6	C8	WHI-P180	23452	RET, KDR
6	C9	NVP-ACC789	23458	VEGFR family
6	C10	RAF709	23820	B-RAF, C-RAF
6	C11	INDY	23895	DYRK1A
6	C12	<i>Unused</i>		
6	D1	<i>Unused</i>		
6	D2	TMB 8 (hydrochloride)	23928	PKC
6	D3	PHT-427	24188	Akt, PDPK1
6	D4	Hesperadin	24199	Multi-kinase
6	D5	Sphingosine (d14:1)	24373	PKC
6	D6	TA-01	24667	CK1 ϵ , CK1 Δ
6	D7	SKI 178	24880	SPHK1, SPHK2
6	D8	XL413 (hydrochloride)	24906	Cdc7
6	D9	Verbascoside	24965	PKC
6	D10	Schizandrin B	24968	ATR
6	D11	ALW-II-41-27	25275	Multi-kinase
6	D12	<i>Unused</i>		
6	E1	<i>Unused</i>		
6	E2	LY3200882	25276	ALK5
6	E3	CP 466,722	25417	ATM
6	E4	A-419259	25605	Src family
6	E5	RK-20448	25606	LCK
6	E6	SGK1 Inhibitor	25652	SGK1, SGK2
6	E7	4-(Dimethylamino)-1-methylpyridinium (iodide)	25861	ChoK
6	E8	Acalisib	25983	PI3K p110 δ
6	E9	LY3177833	25988	Cdc7
6	E10	NIH 12848	26035	PI5P4Ky
6	E11	BMS 794833	26180	c-Met, VEGFR2
6	E12	<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 16 of 30

Plate	Well	Contents	Item Number	Primary Target*
6	F1	Unused		
6	F2	GSK180736A	26182	ROCK-I
6	F3	Pexmetinib	26187	Tie2, p38 α MAPK, p38 β MAPK
6	F4	MK-8745	26190	Aurora A
6	F5	Telatinib	26203	VEGFR family, PDGFR family
6	F6	HMN-214	26210	Plk
6	F7	CH5132799	26410	Class I PI3Ks
6	F8	CMPD101	26808	GRK2, GRK3
6	F9	STK16-IN-1	27370	STK16
6	F10	Streptochlorin	27605	Syk, LYN, Fyn
6	F11	GNE-477	27977	PI3K, mTOR
6	F12	Unused		
6	G1	Unused		
6	G2	D-myo-Inositol-1,3,4-triphosphate (sodium salt)	60972	Ins(3,4,5,6)-P4 kinase
6	G3	APY0201	9001589	PIKfyve
6	G4	A 83-01	9001799	ALK4, ALK5, ALK7
6	G5	PF-299804	9001879	EGFR family
6	G6	WH-4-023	9002067	Src family
6	G7	THZ1	9002215	Cdk7
6	G8	Retreversine	10006606	Negative control
6	G9	Sphingosine Kinase Inhibitor 2	10009222	SPHK1
6	G10	PI3-Kinase α Inhibitor 2	10010177	PI3K p110 α
6	G11	Terreic Acid	10010235	BTK
6	G12	Unused		
6	H1	Unused		
6	H2	AG-825	10010243	HER2
6	H3	AG-17	10010248	EGFR
6	H4	H-8 (hydrochloride)	10010249	PKG
6	H5	RG-14620	10010310	EGFR
6	H6	5-Iodotubercidin	10010375	CK1, PKC, ERK2
6	H7	L-threo-Sphingosine (d18:1)	10010541	PKC
6	H8	PIK-90	10010749	PI3K
6	H9	NH125	10011250	eEF-2K
6	H10	NSC 210902	10011255	CK2
6	H11	CAY10577	10011256	CK2
6	H12	Unused		

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 17 of 30

Plate	Well	Contents	Item Number	Primary Target*
7	A1	Unused		
7	A2	Afatinib	11492	EGFR
7	A3	SL 0101-1	11704	RSK2
7	A4	TBB	14453	CK2
7	A5	DMXAA	14617	VEGFR2
7	A6	Zaltoprofen	14662	PKC
7	A7	Vandetanib	14706	VEGFR2, VEGFR3
7	A8	Ro 3306	15149	Cdk1/cyclin B ₁
7	A9	NU 6027	16177	Cdk1, Cdk2
7	A10	AZD 9291	16237	Mutant EGFR
7	A11	Hemin chloride	16487	eIF2αK1
7	A12	Unused		
7	B1	Unused		
7	B2	CZC-8004	16718	Non-receptor tyrosine kinases
7	B3	CP 724,714	19172	HER2
7	B4	CCT241533 (hydrochloride)	19178	Chk2
7	B5	HG-9-91-01	19179	SIK1, SIK2, SIK3
7	B6	BI-605906	19184	IKKβ
7	B7	PF-3758309	19186	PAK4
7	B8	EW-7197	19231	ALK5
7	B9	Entrectinib	19476	Trk family, ROS1, ALK
7	B10	SU 16f	19555	PDGFRβ
7	B11	Vinpocetine	19578	IKKβ
7	B12	Unused		
7	C1	Unused		
7	C2	HI TOPK 032	19815	TOPK
7	C3	Acalabrutinib	19899	BTK
7	C4	CC-223	19917	mTOR
7	C5	Radotinib	19923	Bcr-Abl
7	C6	CAY10717	19955	Multi-kinase
7	C7	CP21R7	20573	GSK3β
7	C8	DSP-4	20707	Non-kinase inhibitor control
7	C9	FRAX1036	20711	PAK1
7	C10	Dicoumarol	20764	SAPK
7	C11	Daphnetin	20826	EGFR, PKA
7	C12	Unused		

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 18 of 30

Plate	Well	Contents	Item Number	Primary Target*
7	D1	<i>Unused</i>		
7	D2	Necrosulfonamide	20844	MLKL
7	D3	(±)-Necrostatin-2	20924	RIPK1
7	D4	Selonsertib	20972	ASK1
7	D5	eCF309	21118	mTOR
7	D6	DMAT	21182	Pim-1, Pim-3, CK2
7	D7	TNP	21189	IP ₆ K1
7	D8	CDK/CRK Inhibitor	21195	CDKs and CRKs
7	D9	PP121	21196	Multi-kinase
7	D10	R-59-949	21221	DGK-α
7	D11	RSM-932A	21518	CHOKα
7	D12	<i>Unused</i>		
7	E1	<i>Unused</i>		
7	E2	R428	21523	Axl
7	E3	LY2835219	21560	Cdk4, Cdk6
7	E4	SPHINX31	21582	SRPK1
7	E5	AT-13148	21597	Multi-kinase
7	E6	Ro 3280	21669	Plk1
7	E7	βARK1 Inhibitor	21751	GRK2/βARK1
7	E8	AGL 2263	21776	IGF-1R, InsR
7	E9	PI-828	21811	PI3K p110β
7	E10	PKI-166	21896	EGFR
7	E11	MP-A08	21903	SPHK2
7	E12	<i>Unused</i>		
7	F1	<i>Unused</i>		
7	F2	eFT508	21957	MNK1, MNK2
7	F3	YMU1	21981	TMPK
7	F4	XMU-MP-1	22083	MST1, MST2
7	F5	Golvatinib	22143	c-Met, VEGFR2
7	F6	CTX-0294885	22197	Multi-kinase
7	F7	KX1-004	22198	Src
7	F8	LY2090314	22211	GSK3
7	F9	Ki20227	22258	c-FMS
7	F10	N-desethyl Sunitinib	22292	FLK1, PDGFRβ
7	F11	Syringin	22387	AMPKα activator
7	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 19 of 30

Plate	Well	Contents	Item Number	Primary Target*
7	G1	Unused		
7	G2	Nazartinib	22409	Mutant EGFR
7	G3	CC-930	22466	JNK1, JNK2, JNK3
7	G4	BMS 582949	22892	p38α MAPK
7	G5	APY-29	22913	IRE1α
7	G6	R112	22929	Syk
7	G7	ENMD-2076	22956	Multi-kinase
7	G8	WEHI-345	23023	RIPK2
7	G9	Sappanone A	23205	Multi-kinase
7	G10	AG-555	23247	EGFR
7	G11	LY2228820 (mesylate)	23259	p38α MAPK
7	G12	Unused		
7	H1	Unused		
7	H2	SLC5111312 (hydrochloride)	23288	SPHK1, SPHK2
7	H3	SLM6031434	23289	SPHK2
7	H4	GSK872	23300	RIPK3
7	H5	XMD8-87	23354	TNK2
7	H6	Binucleine 2	23367	Aurora B (<i>Drosophila</i>)
7	H7	N ⁶ -Methyladenosine 5'-monophosphate (sodium salt)	23382	Adenylate kinase II
7	H8	Bimiralisib	23441	PI3Kα, mTOR
7	H9	GNE-9605	23446	LRRK2
7	H10	Hexadecyl Methyl Glycerol	60930	PKC
7	H11	Triciribine	10010237	Akt1, Akt2, Akt3
7	H12	Unused		
8	A1	Unused		
8	A2	ABC294640	10587	SPHK2
8	A3	FLLL32	10638	JAK2
8	A4	SC-66	10876	Akt
8	A5	NVP-BKM120	11587	PI3K p110α, PI3K p110β
8	A6	AZD 8330	11588	MEK1, MEK2
8	A7	GDC-0941	11600	PI3Ks
8	A8	SB-505124	11793	ALK5
8	A9	ML-7 (hydrochloride)	11801	MLCK
8	A10	Crizotinib	12087	c-Met, ALK
8	A11	Mubritinib	12096	HER2
8	A12	Unused		

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 20 of 30

Plate	Well	Contents	Item Number	Primary Target*
8	B1	Unused		
8	B2	Tandutinib	12098	PDGFR β , FLT3, c-Kit
8	B3	SU 5402	13182	VEGFR2, FGFR1
8	B4	K252c	13514	PKC
8	B5	NVP-AEW541 (hydrochloride)	13641	IGF-1R
8	B6	AZD 1152 (hydrochloride)	13647	Aurora B
8	B7	CZC-24832	14307	PI3K γ
8	B8	ANA-12	14309	TrkB
8	B9	A-771726	14404	Src family
8	B10	HA-1100 (hydrochloride)	14413	ROCK
8	B11	PF-05212384	14567	PI3K α , PI3K γ , mTOR
8	B12	Unused		
8	C1	Unused		
8	C2	Purvalanol A	14579	Cdc2, Cdk2
8	C3	PF-670462 (hydrochloride)	14588	CK1 ϵ , CK1 δ
8	C4	GSK2578215A	14603	LRRK2, LRRK2 ^{G2019S}
8	C5	TG101209	14696	JAK2, FLT3, RET
8	C6	IPA-3	14759	PAK1
8	C7	SU 9516	14796	Cdk2, Cdk1
8	C8	AS-1892802	14978	ROCKs
8	C9	PF-03814735	15015	Aurora A, Aurora B
8	C10	PF-04708671	15018	S6K1
8	C11	Dovitinib	15220	Multi-kinase
8	C12	Unused		
8	D1	Unused		
8	D2	BI-D1870	15264	RSKs
8	D3	AZ 628	15336	Multi-kinase
8	D4	LY303511 (hydrochloride)	15514	mTOR
8	D5	CCT137690	15552	FLT3, FLT3-ITD, Aurora kinases
8	D6	Akt Inhibitor IV	15569	Akt
8	D7	SL 327	15598	MEK1, MEK2
8	D8	JNK Inhibitor IX	15624	JNK2, JNK3
8	D9	Pelitinib	15627	EGFR
8	D10	ISCK03	15781	c-Kit
8	D11	CID-755673	15924	PKD1, PKD2, PKD3
8	D12	Unused		

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 21 of 30

Plate	Well	Contents	Item Number	Primary Target*
8	E1	Unused		
8	E2	MK2 Inhibitor III	15943	MAPKAPK-2, MAPKAPK-5
8	E3	WHI-P154	16178	EGFR, JAK3
8	E4	AT7519 (hydrochloride)	16231	Cdk2, Cdk4, Cdk5, Cdk9
8	E5	CX-6258	16245	PIMs
8	E6	CYT387	16246	JAK1, JAK2
8	E7	RKI-1447	16278	ROCK-I, ROCK-II
8	E8	TDZD-8	16287	GSK3 β
8	E9	WZ4002	16297	EGFR ^{T790M}
8	E10	XL019	16298	JAK2, JAK2 ^{V617F}
8	E11	XL147 analog 2	16299	PI3K, mTOR
8	E12	Unused		
8	F1	Unused		
8	F2	UNC3230	16401	PIP5K1C
8	F3	SGL-1776	16423	Multi-kinase
8	F4	CHIR124	16553	Chk1
8	F5	LY2784544	16705	JAK2, JAK2 ^{V617F}
8	F6	NVP-BSK805 (hydrochloride)	16714	JAK2, JAK2 ^{V617F}
8	F7	1-Azakenpauillone	16733	GSK3 β
8	F8	R-59-022	16772	DGK
8	F9	CX-4945	16779	CK2, CLK2
8	F10	AT7867	16848	Akt isoforms, S6K1, PKA
8	F11	CFI-400945	16850	Plk4
8	F12	Unused		
8	G1	Unused		
8	G2	N-desmethyl Imatinib	16947	Bcr-Abl
8	G3	ZM 306416	16968	EGFR
8	G4	RAF265	16991	Multi-kinase
8	G5	Binimetinib	16996	MEK1, MEK2
8	G6	BIX02189	16999	MEK5
8	G7	KD 025	17055	ROCK-II
8	G8	BPIQ-I	17224	EGFR
8	G9	AG957	17259	Bcr-Abl, c-Abl
8	G10	IKK2 Inhibitor VI	17276	IKK β
8	G11	GSK2269557	17388	PI3K δ
8	G12	Unused		

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 22 of 30

Plate	Well	Contents	Item Number	Primary Target*
8	H1	<i>Unused</i>		
8	H2	PF-431396	17665	FAK, PYK2
8	H3	KP372-1	17808	Akt
8	H4	AZ 5104	17994	EGFR ^{T790M}
8	H5	Bikinin	18381	ASK kinases
8	H6	BLU-9931	18405	FGFR4
8	H7	AZD 6738	21035	ATR
8	H8	Peficitinib	22954	JAKs
8	H9	THZ531	26386	Cdk12, Cdk13
8	H10	BAY-1895344 (hydrochloride)	26536	ATR
8	H11	KN-92 (hydrochloride)	9000890	Negative control
8	H12	<i>Unused</i>		
9	A1	<i>Unused</i>		
9	A2	Bisindolylmaleimide III	11072	PKC
9	A3	NG 52	17309	Pho85
9	A4	GSK2656157	17372	PERK
9	A5	OSI-027	17379	mTORC1, mTORC2
9	A6	IRAK-1/4 Inhibitor	17540	IRAK1, IRAK4
9	A7	VEGFR2 Kinase Inhibitor II	17544	VEGFR2
9	A8	Adenosine Kinase Inhibitor (hydrate)	17569	Adenosine kinase
9	A9	AZ 20	17589	ATR
9	A10	ML-347	17640	ALK1, ALK2
9	A11	p38 MAPK Inhibitor	17650	p38 MAPK
9	A12	<i>Unused</i>		
9	B1	<i>Unused</i>		
9	B2	GS-9973	17653	Syk
9	B3	VEGFR Tyrosine Kinase Inhibitor II	17654	VEGFR2
9	B4	AMG 706	17671	Multi-kinase
9	B5	JAK2 Inhibitor V	17684	JAK2
9	B6	LH 846	17686	CK1δ
9	B7	AZ191	17693	DYRK1B
9	B8	LDN-212854	17698	ALK2
9	B9	TAK-960	17701	Plk1
9	B10	TTP 22	17702	CK2
9	B11	UNC2250	17703	Mer
9	B12	<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 23 of 30

Plate	Well	Contents	Item Number	Primary Target*
9	C1	<i>Unused</i>		
9	C2	Tie2 Kinase Inhibitor	17858	Tie2
9	C3	PF-477736	17859	Chk1
9	C4	SNS-032	17904	Cdk2
9	C5	3-(4-Pyridyl)indole	17922	ROCK-I
9	C6	Cdk4/6 Inhibitor IV	17974	Cdk4, Cdk6
9	C7	AG-1557	17989	EGFR
9	C8	AVL-292	17993	BTK
9	C9	BS-181 (hydrochloride)	17996	Cdk7
9	C10	Ki8751	18004	VEGFR2
9	C11	XL147	18011	Class I PI3Ks
9	C12	<i>Unused</i>		
9	D1	<i>Unused</i>		
9	D2	ML-167	18098	CLK4
9	D3	GSK461364	18099	Plk1
9	D4	CCT128930	18194	Akt2
9	D5	Ulixertinib (hydrochloride)	18298	ERK1, ERK2
9	D6	A-77-01	18396	ALK5
9	D7	CEP-28122	18406	ALK
9	D8	GDC-0068	18412	Akt1, Akt2, Akt3
9	D9	CGI1746	18414	BTK
9	D10	Rp-8-bromo-Cyclic GMPS (sodium salt)	18441	cGK
9	D11	FLT3 Inhibitor	18462	FLT3
9	D12	<i>Unused</i>		
9	E1	<i>Unused</i>		
9	E2	JNJ-7706621	18494	CDKs, Aurora kinases
9	E3	R788 (sodium salt)	18501	Syk
9	E4	CRT5	18532	PKD1, PKD2, PKD3
9	E5	Akt Inhibitor XI	18604	Akt
9	E6	2-Aminopurine (hydrochloride)	18609	PKR
9	E7	CCT251545	18761	Cdk8, Cdk19
9	E8	4,5,6,7-Tetrabromobenzimidazole	18886	CK2
9	E9	GSK3 β Inhibitor II	19077	GSK3 β
9	E10	LDN-192960 (hydrochloride)	19089	Haspin, DYRK2
9	E11	HTH-01-015	19091	NUAK1
9	E12	<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 24 of 30

Plate	Well	Contents	Item Number	Primary Target*
9	F1	<i>Unused</i>		
9	F2	Purvalanol B	19115	Cdk2/cyclin A, Cdk2/cyclin E, Cdk5/p35, Cdk2/cyclin B
9	F3	TC-S 7001	19129	ROCK-I, ROCK-II
9	F4	MKC-3946	19152	IRE1α
9	F5	CP 673,451	19170	PDGFRα, PDGFRβ
9	F6	PF-4800567	19171	CK1ε
9	F7	GSK269962	19180	ROCK-I, ROCK-II
9	F8	CC-115	19420	mTOR, DNA-PK
9	F9	Narasin (sodium salt)	19447	IκBa phosphorylation
9	F10	YLF-466D	19475	AMPK activator
9	F11	AST-487	19477	Multi-kinase
9	F12	<i>Unused</i>		
9	G1	<i>Unused</i>		
9	G2	RO4987655	19488	MEK
9	G3	MeBIO	19573	Negative control
9	G4	7BIO	19619	FLT3, Aurora C
9	G5	BLZ-945	19626	CSF1R/FMS
9	G6	NPS 1034	19627	Met, Axl
9	G7	Petunidin (chloride)	19755	FAK
9	G8	AP26113	19778	ALK family
9	G9	FIIN-2	19837	FGFR family
9	G10	ON-123300	19902	Multi-kinase
9	G11	Uprosertib	19904	Akt
9	G12	<i>Unused</i>		
9	H1	<i>Unused</i>		
9	H2	GNF-7	19918	Multi-kinase
9	H3	AZD 3759	19931	EGFR
9	H4	HG-10-102-01	20049	LRRK2
9	H5	Cerdulatinib	20076	Syk, JAK family
9	H6	XL388	20237	mTOR
9	H7	LY2603618	20351	Chk1
9	H8	OXA-01	20456	mTORC1, mTORC2
9	H9	PD 168393	21059	EGFR
9	H10	CGP 77675	21089	Src family
9	H11	GDC-0994	21107	ERK1, ERK2
9	H12	<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 25 of 30

Plate	Well	Contents	Item Number	Primary Target*
10	A1	Unused		
10	A2	NU 6140	17271	Cdk2
10	A3	Unused		
10	A4	Tivozanib (hydrate)	18493	VEGFR family, PDGFR family
10	A5	BMS-777607	18517	Met family, TAM family
10	A6	Y11	18521	FAK1
10	A7	Altiratinib	18777	Multi-kinase
10	A8	XMD8-92	19173	ERK5
10	A9	WZ4003	19177	NUAK1, NUAK2
10	A10	AZD 7545	19282	PDK1, PDK2
10	A11	Cobimetinib	19563	MEK1
10	A12	Unused		
10	B1	Unused		
10	B2	Unused		
10	B3	NVS-PAK1-1	19964	PAK1
10	B4	Copanlisib	20354	Class I PI3Ks
10	B5	TX-1918	20840	eEF-2K
10	B6	PKI-179 (hydrochloride)	21202	PI3K, mTOR
10	B7	CYC-116	21205	Aurora A, Aurora B
10	B8	LX7101 (hydrochloride)	21209	LIMKs and ROCKs
10	B9	Fenebrutinib	21346	BTK
10	B10	ZM 323881 (hydrochloride)	21347	VEGFR2
10	B11	NU 2058	21415	Cdk1, Cdk2
10	B12	Unused		
10	C1	Unused		
10	C2	G-5555	21469	Multi-kinase
10	C3	LY2606368	21490	Chk1
10	C4	Cdk1 Inhibitor	21501	Cdk1
10	C5	XL228	21506	Multi-kinase
10	C6	GZD-824	21508	Bcr-Abl
10	C7	Rp-8-bromo-Cyclic AMPS (sodium salt)	21584	PKAs
10	C8	AZD 5438	21598	Multi-kinase
10	C9	Aurora A Inhibitor I	21600	Aurora A
10	C10	BAY 12-17389	21660	Mps1/TTK
10	C11	Fascaplysin (chloride)	21715	Cdk4/cyclin D1
10	C12	Unused		

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 26 of 30

Plate	Well	Contents	Item Number	Primary Target*
10	D1	<i>Unused</i>		
10	D2	SAR407899	21717	ROCKs
10	D3	GSK2292767	21757	PI3Kδ
10	D4	INH6	21772	Nek2
10	D5	GNE-495	21808	MAP4K4
10	D6	JNJ-42756493	21813	FGFR family, VEGFR2
10	D7	Aurora Kinase Inhibitor III	21851	Aurora A
10	D8	Aloisine RP106	21873	Cdk1/cyclinB, Cdk5/p25, GSK3
10	D9	SSR 128129E	22128	FGFR family
10	D10	MPI-0479605	22136	Mps1/TTK
10	D11	G-749	22139	FLT3
10	D12	<i>Unused</i>		
10	E1	<i>Unused</i>		
10	E2	T5601640	22140	LIMK2
10	E3	p38 MAPK Inhibitor IV	22219	p38 MAPKs
10	E4	DB07268	22257	JNK1
10	E5	Pozotinib	22437	EGFR family
10	E6	BLU-554	22438	FGFR4
10	E7	PD 166866	22464	FGFR1
10	E8	AZD 3147	22474	mTOR
10	E9	PRT060318	22476	Syk
10	E10	PF-6274484	22698	EGFR
10	E11	Palomid 529	22706	mTOR
10	E12	<i>Unused</i>		
10	F1	<i>Unused</i>		
10	F2	Epiblastin A	22758	CK1α, CK1ε, CK1δ
10	F3	3-Methylcrotonyl Glycine	22888	Creatine kinase
10	F4	CZC-25146	22931	LRRK2
10	F5	GSK1070916	22953	Aurora B, Aurora C
10	F6	UNC569	22966	TAM family
10	F7	PFK158	22987	PFKFB3
10	F8	FM-381	23204	JAK3
10	F9	5-(Hydroxymethyl)-2'-deoxyuridine	23381	HSV-1 pyrimidine 2'-deoxyribose nucleoside kinase
10	F10	PIK-294	23427	PI3K p110δ
10	F11	CID-797718	23428	PKD1, CAK, Plk1
10	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 27 of 30

Plate	Well	Contents	Item Number	Primary Target*
10	G1	<i>Unused</i>		
10	G2	B-RAF IN 1	23434	B-RAF, C-RAF
10	G3	Solcitinib	23439	JAK1
10	G4	WZ3146	23440	EGFR family
10	G5	Naquotinib	23498	EGFR
10	G6	Brivanib	23690	VEGFR1, VEGFR2, FGFR1
10	G7	TAK-901	23772	Multi-kinase
10	G8	PD 198306	23966	MEK1, MEK2
10	G9	NVP-BAG956	24195	PDK1, Class I PI3Ks
10	G10	AMG 208	24197	c-Met
10	G11	DNA-PK Inhibitor IV	24304	PI3K p110 β , PI3K p110 δ , PI3K p110 γ
10	G12	<i>Unused</i>		
10	H1	<i>Unused</i>		
10	H2	TA-02	24668	CK1 ϵ , CK1 Δ
10	H3	GSK1838705A	24904	IGF-1R, InsR
10	H4	XMD17-109	25448	ERK5
10	H5	CCT241161	25536	Multi-kinase
10	H6	CCT196969	25537	Multi-kinase
10	H7	Compound 43 TAO Kinase Inhibitor	25632	TAOK1, TAOK2
10	H8	CUDC-907	25651	PI3Ks
10	H9	CAY10554	10010301	Cdk5, Cdk2
10	H10	3-O-methyl-N-acetyl-D-Glucosamine	10011056	N-acetylglucosamine kinase, N-acetylmannosamine kinase
10	H11	Hispidin	10012605	PKC β
10	H12	<i>Unused</i>		
11	A1	<i>Unused</i>		
11	A2	K252a	11338	Phosphorylase kinase, PRK1
11	A3	SB 203580	13067	p38 MAPK
11	A4	SU 5416	13342	c-Kit
11	A5	GW 856553X	13614	p38 α MAPK, p38 β MAPK
11	A6	PX 866	13645	PI3K
11	A7	17 β -hydroxy Wortmannin	13812	PI3K
11	A8	(R)-CR8	14006	Cdk2/cyclin E, Cdk2/cyclin A, Cdk1/cyclin B
11	A9	BMS-536924	14666	IGF-1R, InsR
11	A10	Dinaciclib	14707	Cdk1, Cdk2, Cdk5, Cdk9
11	A11	AG-1024	14833	IGF-1R
11	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 28 of 30

Plate	Well	Contents	Item Number	Primary Target*
11	B1	<i>Unused</i>		
11	B2	CID-2011756	15317	PKD
11	B3	MAZ51	16168	VEGFR3
11	B4	Skepinone-L	16974	p38 α MAPK
11	B5	Toyocamycin (hydrate)	17371	RIO1, phosphatidylinositol kinase
11	B6	AMG 319	17383	PI3K δ
11	B7	Cdk4 Inhibitor	17648	Cdk4/cyclin D1
11	B8	BI-847325	18301	Aurora kinases, MEK1, MEK2
11	B9	A-674563	18715	Akt1, PKA, Cdk2
11	B10	GNE-3511	19174	DLK/MAP3K12
11	B11	Mps1-IN-1	19175	Mps1
11	B12	<i>Unused</i>		
11	C1	<i>Unused</i>		
11	C2	Tpl2 Kinase Inhibitor	19710	Tpl2
11	C3	GSK583	19739	RIPK2
11	C4	ETP-46464	19809	ATR, mTOR, DNA-PK
11	C5	MRT67307 (hydrochloride)	19916	Multi-kinase
11	C6	TCS ERK 11e	19932	ERK2
11	C7	INCB 28060	20056	c-Met
11	C8	NPC-15437 (hydrochloride)	20212	PKC
11	C9	BMS-911543	21088	JAK2
11	C10	FLT3 Inhibitor III	21193	FLT3
11	C11	Adenosine 5'-phosphosulfate (sodium salt)	21226	Adenosine 5'-phosphosulfate kinase
11	C12	<i>Unused</i>		
11	D1	<i>Unused</i>		
11	D2	SKLB610	21561	VEGFR2
11	D3	N ⁶ -Benzyladenine	21711	Plant respiratory kinase
11	D4	Compound 401	21769	DNA-PK
11	D5	CFI-402257	21960	Mps1/TTK
11	D6	JH-II-127	22905	LRRK2, LRRK2 ^{G2019S}
11	D7	RKI-1313	22933	ROCK-II
11	D8	GSK2256098	22955	FAK
11	D9	AZD 8931	24196	EGFR, HER2, HER3
11	D10	BI-882370	24273	B-RAF, B-RAFF ^{V600E} , C-RAF (DGF-out)
11	D11	PD 161570	24681	FGFR1
11	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 29 of 30

Plate	Well	Contents	Item Number	Primary Target*
11	E1	Unused		
11	E2	AMG 47a	25339	Multi-kinase
11	E3	Oridonin	25665	Akt1, Akt2
11	E4	CCT244747	25666	Chk1
11	E5	LY2801653	25978	Multi-kinase
11	E6	Flavopiridol	26024	CDKs
11	E7	Unused		
11	E8	Brivanib Alaninate	26081	VEGFR1, VEGFR2, FGFR1
11	E9	TAK-715	26170	p38α MAPK
11	E10	AZD 3463	26172	ALK, IGF-1R
11	E11	GDC-0084	26390	PI3K, mTOR
11	E12	Unused		
11	F1	Unused		
11	F2	IPI-549	26416	PI3Kγ
11	F3	LOXO-101	27056	Trk family
11	F4	Autophinib	27320	VPS34
11	F5	NG 25 (hydrochloride hydrate)	27579	Multi-kinase
11	F6	Hexadecyl Acetyl Glycerol	60920	PKC
11	F7	Jasmonic Acid methyl ester	9000059	Mitochondria-bound hexokinase
11	F8	PI-3065	9002394	PI3K p110δ
11	F9	Genistein	10005167	EGFR, pp50 ^{v-Src} , pp110 ^{bag-fes}
11	F10	D-myo-Inositol-1,3,4,5,6-pentaphosphate (sodium salt)	10007784	PKB/Akt
11	F11	Leelamine (hydrochloride)	10008614	PDK
11	F12	Unused		
11	G1	Unused		
11	G2	CAY10505	10009078	PI3Kγ, CK2
11	G3	PIK-93	10009212	PI4KIIIβ, PI3K p110γ, PI3K p110α
11	G4	SB-415286	10010247	GSK3
11	G5	Lavendustin A	10010268	EGFR
11	G6	AG-213	10010314	EGFR
11	G7	Manumycin A	10010497	IKK
11	G8	HA-1004 (hydrochloride)	10010558	PKG, PKA
11	G9	WHI-P131	10011246	JAK3
11	G10	CAY10576	10011249	IKKε
11	G11	Unused		
11	G12	Unused		

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 30 of 30

Plate	Well	Contents	Item Number	Primary Target*
11	H1	Unused		
11	H2	Unused		
11	H3	Unused		
11	H4	Unused		
11	H5	Unused		
11	H6	Unused		
11	H7	Unused		
11	H8	Unused		
11	H9	Unused		
11	H10	Unused		
11	H11	Unused		
11	H12	Unused		

*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