

# PRODUCT INFORMATION



## Anti-Inflammatory Screening Library

Item No. 31530 • Batch No. 0597693

Page 1 of 10

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

Plate	Well	Contents	Item Number	Primary Target*
1	A1	Unused		
1	A2	10-Nitrolinoleate	10037	PPAR $\gamma$
1	A3	FR122047 (hydrochloride)	10039	COX-1
1	A4	BM 567	10155	TXA $_2$ Synthase, TP Receptor
1	A5	Doramapimod	10460	p38 MAPK
1	A6	CAY10650	10743	cPLA $_2\alpha$
1	A7	CAY10649	10804	5-LO
1	A8	HZ52	10888	5-LO
1	A9	Dexamethasone	11015	Anti-inflammatory Activity/ Glucocorticoid
1	A10	TPPU	11120	sEH
1	A11	R406	11422	Syk
1	A12	Unused		
1	B1	Unused		
1	B2	Ruxolitinib	11609	JAK1, JAK2
1	B3	1-Naphthyl 3,5-dinitrobenzoate	11729	5-LO, mPGES-1
1	B4	Treprostinil (diethanolamine salt)	11927	IP $_2$ , DP $_2$ , and EP $_2$ Receptors
1	B5	Fingolimod	11975	Immunomodulatory Activity
1	B6	Cyclosporin A	12088	Immunosuppressant/Calcineurin
1	B7	PD 0325901	13034	MEK
1	B8	Tranilast	13044	Anti-inflammatory and Immunomodulatory Activity
1	B9	CAY10589	13164	5-LO, mPGES-1
1	B10	CP 80,633	13183	PDE4
1	B11	MF63	13217	mPGES-1
1	B12	Unused		
1	C1	Unused		
1	C2	ZLJ-6	13271	COX-1, COX-2, 5-LO
1	C3	Pyrophenone	13294	cPLA $_2\alpha$
1	C4	CGP 57380	13322	MNK1
1	C5	PPM-18	13327	NF- $\kappa$ B Signaling
1	C6	CAY10606	13381	5-LO
1	C7	GW 856553X	13614	p38 $\alpha$ MAPK, p38 $\beta$ MAPK
1	C8	CAY10614	13615	TLR4
1	C9	YS-121	13665	5-LO, mPGES-1
1	C10	GW 1929	13689	PPAR $\gamma$
1	C11	AX 048	13823	cPLA $_2$
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, 03/18/2022

#### 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 10

Plate	Well	Contents	Item Number	Primary Target*
1	D1	<i>Unused</i>		
1	D2	TAK-242	13871	TLR4
1	D3	Methotrexate (hydrate)	13960	Anti-inflammatory Activity/ Nucleic Acid Signaling
1	D4	Azathioprine	13987	Immunosuppressant/ Nucleic Acid Signaling
1	D5	S18886	14059	TP Receptor
1	D6	Amlexanox	14181	Anti-inflammatory Activity, IKK $\epsilon$ , TBK1
1	D7	Chloroquine (phosphate)	14194	Autophagy, Antimalarial
1	D8	Tanshinone IIA	14323	General Anti-inflammatory Activity
1	D9	S3I-201	14336	STAT3
1	D10	MK2 Inhibitor IV	14399	MK2
1	D11	Melatonin	14427	Antioxidant Activity
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	SCH 79797 (hydrochloride)	14633	PAR1
1	E3	Ginkgolide B	14636	PAFR
1	E4	L-Ascorbic Acid	14656	Antioxidant Activity
1	E5	WP1066	14736	STAT3
1	E6	Leflunomide	14860	Prodrug, T Cell Activation/ Pyrimidine Synthesis
1	E7	Meloxicam	14906	COX-2
1	E8	BX-795	14932	Multi-kinase
1	E9	Azithromycin	15004	Antibiotic, Anti-inflammatory Activity
1	E10	Methylprednisolone	15013	Immunosuppressant/ Glucocorticoid
1	E11	Prazosin (hydrochloride)	15023	$\alpha_1$ -Adrenergic Receptors
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	JSH-23	15036	NF- $\kappa$ B Signaling
1	F3	TPCA-1	15115	IKK $\beta$
1	F4	CAY10678	15129	mPGES-1
1	F5	LCK Inhibitor	15135	LCK
1	F6	CU CPT 22	15244	TLR1/TLR2 Heterodimers
1	F7	Auranofin	15316	Anti-inflammatory Activity, Immunosuppressant
1	F8	Ebastine	15372	Histamine H <sub>1</sub> Receptor
1	F9	Ro 106-9920	15373	NF- $\kappa$ B Signaling
1	F10	Budesonide	15407	Anti-inflammatory Activity/ Glucocorticoid
1	F11	Ticagrelor	15425	Purinergic P2Y <sub>12</sub> Receptor
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 10

Plate	Well	Contents	Item Number	Primary Target*
1	G1	<i>Unused</i>		
1	G2	Pirinixic Acid Aminothiazole	15483	5-LO, mPGES-1
1	G3	BW 755C	15490	COX-1, COX-2, 5-LO
1	G4	FR180204	15544	ERK1, ERK2
1	G5	CAY10682	15574	NF- $\kappa$ B Signaling
1	G6	MK2 Inhibitor III	15943	MK2
1	G7	5,15-DPP	16090	STAT3
1	G8	ATB-346	16234	Anti-inflammatory Activity/ NSAID
1	G9	Ibrutinib	16274	BTK
1	G10	TG101348	16289	JAK2, FLT3, RET
1	G11	Trametinib	16292	MEK1, MEK2
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	trans-AUCB	16568	sEH
1	H3	BMS 345541 (trifluoroacetate salt)	16667	IKK $\alpha$ , IKK $\beta$
1	H4	Baricitinib	16707	JAK1, JAK2
1	H5	R-7050	16870	TNF- $\alpha$ Receptor Signaling
1	H6	Binimetinib	16996	MEK1, MEK2
1	H7	TAK-733	16998	MEK1
1	H8	BIX02189	16999	MEK5, ERK5
1	H9	BB-Cl-Amidine	17079	NETosis/PAD
1	H10	CGS 21680 (hydrochloride)	17126	Adenosine A <sub>2A</sub> Receptor
1	H11	IMD 0354	17290	IKK $\beta$
1	H12	<i>Unused</i>		
2	A1	<i>Unused</i>		
2	A2	Antileukinate	17359	CXCR2
2	A3	Tenidap	17413	COX-1
2	A4	(5Z)-7-Oxozeaenol	17459	TAK1
2	A5	GSK484 (hydrochloride)	17488	NETosis/PAD4
2	A6	TNF- $\alpha$ Antagonist	17492	TNF- $\alpha$ Signaling
2	A7	NF- $\kappa$ B Inhibitor (trifluoroacetate salt)	17493	NF- $\kappa$ B Signaling
2	A8	MCC950	17510	NLRP3
2	A9	IRAK-1/4 Inhibitor	17540	IRAK1, IRAK4
2	A10	BIIB-057	17652	Syk
2	A11	Sivelestat (sodium salt hydrate)	17779	NETosis/Neutrophil Elastase
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 10

Plate	Well	Contents	Item Number	Primary Target*
2	B1	<i>Unused</i>		
2	B2	Hydroxychloroquine (sulfate)	17911	Autophagy, Antimalarial
2	B3	JNK Inhibitor XVI	18096	JNK1, JNK2, JNK3
2	B4	Dipyridamole	18189	PDE5A
2	B5	Fexofenadine (hydrochloride)	18191	Histamine H <sub>1</sub> Receptor
2	B6	Syk Inhibitor II	18202	Syk
2	B7	Ciclesonide	18227	Anti-inflammatory Activity/ Glucocorticoid Prodrug
2	B8	Piriprost (potassium salt)	18228	5-LO
2	B9	BX 471	18503	CCR1
2	B10	TAPI-1	18505	ADAM17/TACE
2	B11	TLR4-C34	18512	TLR4 Signaling
2	B12	<i>Unused</i>		
2	C1	<i>Unused</i>		
2	C2	BAY-678	18615	NETosis/Neutrophil Elastase
2	C3	Oclacitinib	18722	JAK Family
2	C4	CAY10700	18905	mPGES-1
2	C5	BMX-IN-1	19165	BMX, BTK
2	C6	SCH 772984	19166	ERK1, ERK2
2	C7	Arglabin	19502	Anti-inflammatory Activity/ NLRP3 Inflammasome
2	C8	Tpl2 Kinase Inhibitor	19710	Tpl2
2	C9	Acalabrutinib	19899	BTK
2	C10	MRT67307 (hydrochloride)	19916	Multi-kinase
2	C11	Ozanimod	19922	S1P <sub>1</sub> and S1P <sub>5</sub>
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	N-acetyl-L-Cysteine	20261	Antioxidant Activity
2	D3	Fluticasone Propionate	20703	Anti-inflammatory Activity/ Corticosteroid
2	D4	Hydrocortisone	20739	Anti-inflammatory Activity/ Glucocorticoid
2	D5	(±)-Necrostatin-2	20924	RIPK1
2	D6	Danirixin	20980	CXCR2
2	D7	(S)-(+)-Clopidogrel (sulfate)	21002	Prodrug, Purinergic P2Y <sub>12</sub> Receptor
2	D8	BAY 59-3074	21036	CB <sub>1</sub> and CB <sub>2</sub> Receptors
2	D9	IL-1R Antagonist	21349	IL-1RI
2	D10	Mometasone Furoate	21365	Anti-inflammatory Activity/ Glucocorticoid Prodrug
2	D11	Bentamapimod	21482	JNK1, JNK2, JNK3
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 10

Plate	Well	Contents	Item Number	Primary Target*
2	E1	<i>Unused</i>		
2	E2	R428	21523	Axl
2	E3	Mycophenolic Acid	21716	Immunosuppressant/IMPDH
2	E4	MK-886	21753	FLAP
2	E5	SC-75741	21818	NF-κB Signaling
2	E6	eFT508	21957	MNK1, MNK2
2	E7	PF-4136309	21965	CCR2
2	E8	Ponesimod	22053	S1P <sub>1</sub>
2	E9	PF-1355	22222	Myeloperoxidase
2	E10	Timosaponin AIII	22236	General Anti-Inflammatory Activity
2	E11	Clonixin	22284	Anti-inflammatory/NSAID
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	IT-901	22298	NF-κB Signaling
2	F3	CC-930	22466	JNK1, JNK2, JNK3
2	F4	R-8507	22479	TNF-αRI
2	F5	PF-4693627	22768	mPGES-1
2	F6	T-5224	22904	c-Fos/AP-1
2	F7	WEHI-345	23023	RIPK2
2	F8	Fluocinolone acetonide	23027	Anti-inflammatory Activity/ Corticosteroid
2	F9	Vorapaxar	23119	PAR1
2	F10	Mizoribine	23128	Immunosuppressant
2	F11	LY2228820 (mesylate)	23259	p38α MAPK
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	GSK872	23300	RIPK3
2	G3	Delanzomib	23557	Cymotrypsin-like Proteasome, NF-κB Signaling
2	G4	Halcinonide	23780	Anti-inflammatory Activity/ Glucocorticoid
2	G5	Flunisolide	23842	Anti-inflammatory Activity/ Corticosteroid
2	G6	Ceralifimod	24159	S1P <sub>1</sub> and S1P <sub>5</sub>
2	G7	Takinib	24161	TAK1
2	G8	Flumethasone	24184	Anti-inflammatory Activity/ Corticosteroid
2	G9	Dapansutrile	24671	NLRP3 Inflammasome
2	G10	Cryogenine	25287	Prostaglandin Synthetase
2	G11	AMG 47a	25339	Multi-kinase
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 10

Plate	Well	Contents	Item Number	Primary Target*
2	H1	<i>Unused</i>		
2	H2	CU-CPT8m	25349	TLR8
2	H3	Acumapimod	25832	p38α MAPK
2	H4	(±)-α-Tocopherol	25985	Antioxidant Activity
2	H5	AZD 9668	26083	NETosis/Neutrophil Elastase
2	H6	GS-143	26138	NF-κB Signaling
2	H7	Verdinexor	26171	XPO1, Antiviral
2	H8	Lipoxygenin	26310	5-LO
2	H9	TAK-779	26744	CCR2b, CCR5, CXCR3
2	H10	Linarin	26900	General Anti-inflammatory Activity
2	H11	(±)-γ-Tocopherol	27193	Antioxidant and Anti-inflammatory Activity
2	H12	<i>Unused</i>		
3	A1	<i>Unused</i>		
3	A2	Oroxlylin A	27363	General Anti-inflammatory Activity
3	A3	NG 25 (hydrochloride hydrate)	27579	Multi-kinase
3	A4	CA-4948	27598	IRAK4
3	A5	Diflapolin	28103	FLAP, sEH
3	A6	Fraxin	28217	General Anti-inflammatory Activity
3	A7	GI 254023X	28284	ADAM10
3	A8	AZD 5069	28297	CXCR2
3	A9	BP-1-102	28368	STAT3
3	A10	<i>Unused</i>		
3	A11	INCB 3344	28433	CCR2
3	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 7 of 10

Plate	Well	Contents	Item Number	Primary Target*
3	B1	<i>Unused</i>		
3	B2	NBI 74330	28442	CXCR3
3	B3	INF39	28476	NLRP3 Inflammasome
3	B4	CU-CPT9a	28722	TLR8
3	B5	6(5H)-Phenanthridinone	28730	PARP1, PARP2
3	B6	TP-0903	28757	Axl, Aurora A, Aurora B
3	B7	IM-93	28794	NETosis
3	B8	Tenuigenin	28824	General Anti-inflammatory Activity
3	B9	SRI-011381	28914	TGF- $\beta$ Signaling
3	B10	Zanubrutinib	28924	BTK
3	B11	Evobrutinib	29029	BTK
3	B12	<i>Unused</i>		
3	C1	<i>Unused</i>		
3	C2	$\alpha$ -Mangostin	29108	General Anti-inflammatory Activity
3	C3	SNS-062 analog	29200	BTK
3	C4	GSK2982772	29230	RIPK1
3	C5	Wogonoside	29332	General Anti-inflammatory Activity
3	C6	Merimepodib	29351	Immunosuppressant/IMPDH
3	C7	Prunetin	29432	General Anti-inflammatory Activity
3	C8	LDC-1267	29459	Axl, Mer, Tyro3
3	C9	YM-976	29529	PDE4
3	C10	Dehydrolithocholic Acid	29544	TGR5, VDR, FXR, PXR, ROR $\gamma$ t
3	C11	Senicapoc	29679	IK $_{Ca}$ 1/K $_{Ca}$ 3.1 Channels
3	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 8 of 10

Plate	Well	Contents	Item Number	Primary Target*
3	D1	<i>Unused</i>		
3	D2	SC144 (hydrochloride)	29714	gp130
3	D3	Betamethasone 21-phosphate (sodium salt hydrate)	29778	Anti-inflammatory Activity/ Glucocorticoid
3	D4	CU-115	29815	TLR8
3	D5	Fluticasone Furoate	29878	Anti-inflammatory Activity/ Glucocorticoid
3	D6	ML-209	30164	ROR $\gamma$ t
3	D7	TG11-77 (hydrochloride)	30188	EP $_2$ Receptor
3	D8	Lapachol	30190	General Anti-inflammatory Activity
3	D9	NLRP3i	30895	NLRP3 Inflammasome
3	D10	CU-CPT4a	30951	TLR3
3	D11	Parthenolide	70080	NF- $\kappa$ B Signaling
3	D12	<i>Unused</i>		
3	E1	<i>Unused</i>		
3	E2	Indomethacin	70270	COX-1, COX-2
3	E3	( $\pm$ )-Ibuprofen	70280	COX-1, COX-2
3	E4	(+)-Naproxen	70290	COX-1, COX-2
3	E5	SC-560	70340	COX-1
3	E6	Tolfenamic Acid	70480	COX-2
3	E7	NS-398	70590	COX-2
3	E8	Caffeic Acid	70602	5-LO, 12-LO
3	E9	CAY10416	70635	COX-2, 5-LO
3	E10	DuP-697	70645	COX-2
3	E11	<i>trans</i> -Resveratrol	70675	Antioxidant and Anti-inflammatory Activity
3	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 9 of 10

Plate	Well	Contents	Item Number	Primary Target*
3	F1	<i>Unused</i>		
3	F2	Caffeic Acid phenylethyl ester	70750	NF-κB Signaling
3	F3	Rosiglitazone	71740	PPARγ
3	F4	Curcumin	81025	Antioxidant and Anti-inflammatory Activity
3	F5	Colchicine	9000760	Microtubule Polymerization, Anti-Inflammatory Activity
3	F6	Ketorolac	9001148	COX-1, COX-2
3	F7	APY0201	9001589	PIKFYVE
3	F8	AF3485	9001962	mPGES-1
3	F9	WH-4-023	9002067	LCK, Src
3	F10	Luteolin	10004161	General Anti-inflammatory Activity, 15-LO
3	F11	SC-236	10004219	COX-2
3	F12	<i>Unused</i>		
3	G1	<i>Unused</i>		
3	G2	cis-Resveratrol	10004235	Antioxidant and Anti-inflammatory Activity
3	G3	N,N'-Dicyclohexylurea	10004971	sEH
3	G4	Quercetin (hydrate)	10005169	Antioxidant and Anti-inflammatory Activity
3	G5	QNZ	10006734	NF-κB Signaling
3	G6	Zileuton	10006967	5-LO
3	G7	Licofelone	10007692	COX, 5-LO
3	G8	CUDA	10007923	sEH
3	G9	FK-506	10007965	Immunosuppressant
3	G10	SPD-304	10008012	TNF-α Signaling
3	G11	α-Tocotrienol	10008377	Antioxidant Activity
3	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 10 of 10

Plate	Well	Contents	Item Number	Primary Target*
3	H1	<i>Unused</i>		
3	H2	Rofecoxib	10010260	COX-2
3	H3	GW 842166X	10010372	CB <sub>2</sub> Receptor
3	H4	CJ-42794	10010428	EP <sub>4</sub> Receptor
3	H5	SP 600125	10010466	JNK1, JNK2, JNK3
3	H6	CAY10576	10011249	IKKε
3	H7	<i>Unused</i>		
3	H7	<i>Unused</i>		
3	H9	<i>Unused</i>		
3	H10	<i>Unused</i>		
3	H11	<i>Unused</i>		
3	H12	<i>Unused</i>		

\*Primary Target is based on available IC<sub>50</sub> and/or K<sub>i</sub> 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