

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 γ
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- κ B Signaling
1	C6	CAY10606	13381	5-LO
1	C7	GW 856553X	13614	p38 α MAPK, p38 β MAPK
1	C8	CAY10614	13615	TLR4
1	C9	YS-121	13665	5-LO, mPGES-1
1	C10	GW 1929	13689	PPAR γ
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, 09/04/2024

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