

PRODUCT INFORMATION



Pesticide Screening Library

Item No. 36390 • Batch No. 0649880

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

Page 1 of 7

Plate	Well	Contents	Item Number	Activity
1	A1	Unused		
1	A2	2,4-Dichlorophenoxy Acetic Acid	10190	Herbicide
1	A3	Pellitorine	11662	Insecticide
1	A4	Ethyl <i>p</i> -methoxycinnamate	11740	Insecticide
1	A5	Asperlactone	14435	Insecticide
1	A6	Rocaglamide	14841	Insecticide
1	A7	Validamycin A	15326	Fungicide
1	A8	Venturicidin A	15377	Fungicide
1	A9	Thaxtomin A	15601	Herbicide
1	A10	2,4-Diacetylphloroglucinol	16345	Fungicide
1	A11	Spinosyn A	16528	Insecticide
1	A12	Unused		
1	B1	Unused		
1	B2	Unused		
1	B3	(S)-(+)-Methoprene	16807	Insecticide
1	B4	Diquat-d ₄ (bromide)	18250	Herbicide
1	B5	Milbemycin A ₃	18360	Acaricide, Insecticide
1	B6	Propiconazole	18853	Fungicide
1	B7	Paclobutrazol	18864	Pesticide
1	B8	Unused		
1	B9	Pyoluteorin	19457	Herbicide
1	B10	Cyclosporin B	19525	Insecticide
1	B11	7-methoxy Apigeninidin (chloride)	19757	Fungicide
1	B12	Unused		
1	C1	Unused		
1	C2	6',7'-Epoxybergamottin	19760	Insecticide
1	C3	Sakuranetin	19761	Fungicide
1	C4	Simetryn	20392	Herbicide
1	C5	Fenpropimorph	20875	Fungicide
1	C6	Nitenpyram	21271	Insecticide
1	C7	Manzamine A	21444	Insecticide
1	C8	Fluensulfone	21554	Nematicide
1	C9	α -Humulene	21574	Insecticide
1	C10	α -Pinene	21576	Insecticide
1	C11	(-)- β -Pinene	21577	Insecticide
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, 06/14/2023

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 7

Plate	Well	Contents	Item Number	Activity
1	D1	<i>Unused</i>		
1	D2	Dianemycin	21679	Insecticide
1	D3	Pyriproxyfen	21872	Insecticide
1	D4	Azadirachtin	21893	Insecticide
1	D5	(±)-Dunnione	21943	Fungicide
1	D6	Avermectin B _{1a}	22000	Insecticide
1	D7	Fluralaner	22061	Acaricide
1	D8	Adipostatin A	22163	Insecticide
1	D9	Fluazuron	22283	Acaricide
1	D10	<i>Unused</i>		
1	D11	Diflubenzuron	23095	Insecticide
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	(-)-Guaiol	23172	Insecticide
1	E3	(1R)-(+)-Camphor	23175	Pesticide
1	E4	(1S)-(-)-Camphor	23176	Acaricide
1	E5	α-Terpinene	23178	Acaricide
1	E6	Tetrachlorohydroquinone	23231	Pesticide
1	E7	Isoxanthohumol	23233	Pesticide
1	E8	Amiprofos-methyl	23379	Herbicide
1	E9	Methazolamide	23413	Insecticide
1	E10	Herbicidin A	23485	Herbicide
1	E11	Brevianamide F	23493	Pesticide
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	Cellocidin	23637	Pesticide
1	F3	Nornidulin	23646	Insecticide
1	F4	Acephate	24038	Insecticide
1	F5	Diuron	24040	Herbicide
1	F6	Prochloraz	24051	Pesticide
1	F7	Bitertanol	24054	Fungicide
1	F8	Cypermethrin	24117	Insecticide
1	F9	Hexaconazole	24118	Fungicide
1	F10	Fipronil	24128	Pesticide
1	F11	Chlorantraniliprole	24140	Insecticide
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 7

Plate	Well	Contents	Item Number	Activity
1	G1	<i>Unused</i>		
1	G2	Chlorfenapyr	24141	Insecticide
1	G3	Chlorothalonil	24142	Fungicide
1	G4	3-hydroxy Carbofuran	24146	Insecticide
1	G5	Aldicarb sulfone	24210	Insecticide
1	G6	Aldicarb Sulfoxide	24211	Insecticide
1	G7	Acibenzolar-S-methyl	24212	Fungicide
1	G8	Chlorotoluron	24225	Herbicide
1	G9	Cyfluthrin	24231	Insecticide
1	G10	<i>o,p'</i> -DDE	24235	Pesticide
1	G11	<i>p,p'</i> -DDE	24241	Pesticide
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	<i>o,p'</i> -DDT	24242	Insecticide
1	H3	Diniconazole	24248	Fungicide
1	H4	Endosulfan I	24253	Insecticide
1	H5	Endosulfan II	24254	Insecticide
1	H6	Endosulfan sulfate	24255	Insecticide
1	H7	(±)-Camphene	24478	Pesticide
1	H8	(+)-Nootkatone	24910	Insecticide
1	H9	Cyclo(L-Leu-L-Pro)	24941	Pesticide
1	H10	Herboxidiene	25136	Herbicide
1	H11	Fenoxycarb	25607	Insecticide
1	H12	<i>Unused</i>		
2	A1	<i>Unused</i>		
2	A2	Flonicamid	25608	Insecticide
2	A3	(E)-Fenpyroximate	25625	Acaricide
2	A4	Methiocarb	25627	Acaricide, Insecticide
2	A5	Propoxur	25628	Insecticide
2	A6	Spirotetramat	25634	Insecticide
2	A7	<i>Unused</i>		
2	A8	<i>trans</i> -Nerolidol	25778	Insecticide
2	A9	Etofenprox	25785	Insecticide
2	A10	Phosmet	25812	Acaricide, Insecticide
2	A11	Kresoxim-methyl	25816	Fungicide
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 7

Plate	Well	Contents	Item Number	Activity
2	B1	<i>Unused</i>		
2	B2	Spiromesifen	25822	Acaricide, Insecticide
2	B3	2-chloro Stearic Acid	26108	Pesticide
2	B4	Anacardic Acid triene	26611	Pesticide
2	B5	Spinetoram	26623	Insecticide
2	B6	Undecanoic Acid methyl ester	26720	Insecticide
2	B7	Bromadiolone	26854	Rodenticide
2	B8	Cyantraniliprole	27309	Insecticide
2	B9	<i>Unused</i>		
2	B10	Aurantioobtusin	27481	Insecticide
2	B11	NG 012	27509	Fungicide
2	B12	<i>Unused</i>		
2	C1	<i>Unused</i>		
2	C2	OM173-αA	27570	Fungicide
2	C3	Ethoprop	27609	Insecticide
2	C4	Ginkgoneolic Acid	27656	Pesticide
2	C5	Virustomycin A	27736	Fungicide
2	C6	<i>Unused</i>		
2	C7	Methiocarb-d ₃	27776	Fungicide
2	C8	Pogostone	27812	Insecticide
2	C9	Cymoxanil-d ₃	27840	Fungicide
2	C10	Umbelliprenin	27894	Insecticide
2	C11	Schaftoside	27950	Pesticide
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	Averantin	28059	Fungicide
2	D3	1-Hydroxyphenazine	28064	Fungicide
2	D4	Eprinomectin	28182	Acaricide, Insecticide
2	D5	Precocene II	28480	Insecticide
2	D6	Phomalactone	28854	Fungicide
2	D7	<i>Unused</i>		
2	D8	5,7-Dihydroxychromone	29006	Fungicide
2	D9	(±)-Indoxacarb	29165	Insecticide
2	D10	Oosporein	29173	Fungicide
2	D11	Fenvalerate	29399	Acaricide, Insecticide
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 7

Plate	Well	Contents	Item Number	Activity
2	E1	<i>Unused</i>		
2	E2	Monocerin	29572	Fungicide, Insecticide
2	E3	Pyrenocine A	29622	Fungicide
2	E4	Tropolone	29674	Fungicide
2	E5	Estragole	29824	Insecticide
2	E6	(S)-(-)-Perillaldehyde	29872	Pesticide
2	E7	Nimbin	29884	Fungicide, Insecticide
2	E8	(4'-Hydroxy)phenoxybenzoic Acid	29891	Insecticide
2	E9	2-Hydroxy-4-methoxybenzaldehyde	30091	Fungicide
2	E10	(-)-Menthol	30384	Pesticide
2	E11	Abametapir	30684	Insecticide
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	<i>Unused</i>		
2	F3	Triflumuron	30858	Insecticide
2	F4	Mahanimbine	30897	Insecticide
2	F5	Chlorpyrifos-d ₁₀	30953	Insecticide
2	F6	Dehydrocorydaline (chloride)	30972	Herbicide
2	F7	Dihydrosanguinarine	31139	Fungicide
2	F8	Anagyrine	31169	Pesticide
2	F9	Fumitremorgin B	31173	Fungicide, Insecticide
2	F10	Propiconazole-d ₇	31177	Fungicide
2	F11	Sclareolide	31371	Fungicide
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	Ascochin	31450	Fungicide
2	G3	Alyssin	31513	Fungicide
2	G4	Diocanoyl Glycol	31515	Insecticide
2	G5	TFM	32870	Piscicide
2	G6	3-Oxobetulin	33066	Fungicide
2	G7	Tebuconazole-d ₉	33287	Fungicide
2	G8	Afoxolaner	33317	Acaricide, Insecticide
2	G9	7,10-dihydroxy-8(E)-Octadecenoic Acid	33447	Pesticide
2	G10	Piperitone	33490	Insecticide
2	G11	N,N'-Diphenethylurea	33580	Fungicide
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 7

Plate	Well	Contents	Item Number	Activity
2	H1	<i>Unused</i>		
2	H2	Tubercycin B	33659	Fungicide
2	H3	Octahydrocurcumin	33686	Fungicide
2	H4	Piericidin B	33721	Insecticide
2	H5	Dinotefuran	33802	Insecticide
2	H6	(S)-Canadine	33805	Insecticide
2	H7	Carveol	33831	Pesticide
2	H8	Methazolamide-d ₆	33933	Insecticide
2	H9	Luteolin 7-O-Rutinoside	34035	Fungicide
2	H10	Toosendanin	34042	Insecticide
2	H11	Isoscopoletin	34117	Fungicide
2	H12	<i>Unused</i>		
3	A1	<i>Unused</i>		
3	A2	Yunaconitine	34127	Insecticide
3	A3	Terbuthylazine	34192	Herbicide
3	A4	Loliolide	34196	Pesticide
3	A5	Naptalam	34285	Herbicide
3	A6	Quassin	34382	Insecticide
3	A7	(+)- δ -Cadinene	34444	Insecticide
3	A8	Glufosinate-d ₃ (hydrochloride)	34447	Herbicide
3	A9	Tuberostemonine	34457	Insecticide
3	A10	Dichlorvos-d ₆	34517	Insecticide
3	A11	Pendimethalin	34557	Herbicide
3	A12	<i>Unused</i>		
3	B1	<i>Unused</i>		
3	B2	Arecaidine propargyl ester (hydrobromide)	34568	Insecticide
3	B3	Menthone	34683	Insecticide
3	B4	4-Acetamidobenzenesulfonamide	34689	Herbicide
3	B5	Allocryptopine	34751	Fungicide
3	B6	N-Methylcytisine	34752	Pesticide
3	B7	Murrayone	34767	Fungicide
3	B8	Simazine-d ₁₀	35209	Herbicide
3	B9	Nanaomycin A	9001479	Fungicide
3	B10	Vinclozolin M ₂	10007452	Fungicide
3	B11	(+)-Goniothalesdiol	10008886	Pesticide
3	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 7 of 7

Plate	Well	Contents	Item Number	Activity
3	C1	<i>Unused</i>		
3	C2	(-)-Deguelin	10010706	Insecticide
3	C3	4-phenyl-1,2,3-Thiadiazole	10011055	Fungicide, Herbicide
3	C4	<i>Unused</i>		
3	C5	<i>Unused</i>		
3	C6	<i>Unused</i>		
3	C7	<i>Unused</i>		
3	C8	<i>Unused</i>		
3	C9	<i>Unused</i>		
3	C10	<i>Unused</i>		
3	C11	<i>Unused</i>		
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