

PRODUCT INFORMATION



Microbiome Metabolite Screening Library

Item No. 36305 • Batch No. 0638687

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

Page 1 of 8

Plate	Well	Contents	Item Number	Class
1	A1	Unused		
1	A2	Trimethylamine N-oxide	17354	Amine
1	A3	Acetaminophen	10024	Aromatic amine
1	A4	Unused		
1	A5	Unused		
1	A6	Unused		
1	A7	Unused		
1	A8	Unused		
1	A9	Unused		
1	A10	Unused		
1	A11	Unused		
1	A12	Unused		
1	B1	Unused		
1	B2	Spermidine	14918	Amino acid derivative
1	B3	Spermine	18041	Amino acid derivative
1	B4	Unused		
1	B5	Unused		
1	B6	Betaine	26814	Amino acid derivative
1	B7	(±)-Synephrine (hydrochloride)	28398	Amino acid derivative
1	B8	Unused		
1	B9	Unused		
1	B10	N,N-Dimethylglycine (hydrochloride)	33830	Amino acid derivative
1	B11	N-Acetyl-DL-Aspartic Acid	34635	Amino acid derivative
1	B12	Unused		
1	C1	Unused		
1	C2	Glycoyamine	35792	Amino acid derivative
1	C3	N ^G ,N ^G -dimethyl-L-Arginine (hydrochloride)	80230	Amino acid derivative
1	C4	Unused		
1	C5	Unused		
1	C6	Unused		
1	C7	Unused		
1	C8	Unused		
1	C9	Unused		
1	C10	Unused		
1	C11	Unused		
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, 12/14/2023

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 8

Plate	Well	Contents	Item Number	Class
1	D1	<i>Unused</i>		
1	D2	L-Kynurenine	11305	Aromatic amino acid derivative
1	D3	5-Methoxyindole-3-acetic acid	14019	Aromatic amino acid derivative
1	D4	(±)- <i>para</i> -Octopamine (hydrochloride)	14279	Aromatic amino acid derivative
1	D5	Serotonin (hydrochloride)	14332	Aromatic amino acid derivative
1	D6	Metanephrine	15408	Aromatic amino acid derivative
1	D7	<i>trans</i> -Urocanic Acid	16228	Aromatic amino acid derivative
1	D8	Phenylacetyl L-Glutamine	16724	Aromatic amino acid derivative
1	D9	Kynurenic Acid (hydrate)	16792	Aromatic amino acid derivative
1	D10	Indoxyl Sulfate (potassium salt)	16926	Aromatic amino acid derivative
1	D11	Tyramine	18601	Aromatic amino acid derivative
1	D12	<i>Unused</i>		
1	E1	<i>Unused</i>		
1	E2	Phenylacetic Acid	18709	Aromatic amino acid derivative
1	E3	Xanthurenic Acid	19191	Aromatic amino acid derivative
1	E4	3-hydroxy Anthranilic Acid	20512	Aromatic amino acid derivative
1	E5	<i>p</i> -Coumaric Acid	20929	Aromatic amino acid derivative
1	E6	Tryptamine	20995	Aromatic amino acid derivative
1	E7	Dopamine (hydrochloride)	21992	Aromatic amino acid derivative
1	E8	5-hydroxy Indole-3-acetic Acid	22889	Aromatic amino acid derivative
1	E9	4-Aminohippuric Acid	23726	Aromatic amino acid derivative
1	E10	Oxindole	23851	Aromatic amino acid derivative
1	E11	3,4-Dihydroxyphenylacetic Acid	24912	Aromatic amino acid derivative
1	E12	<i>Unused</i>		
1	F1	<i>Unused</i>		
1	F2	6-OHDA (hydrobromide)	25330	Aromatic amino acid derivative
1	F3	4-Hydroxyphenylacetic Acid	26613	Aromatic amino acid derivative
1	F4	Dihydrocaffeic Acid	28390	Aromatic amino acid derivative
1	F5	Tryptophol	28404	Aromatic amino acid derivative
1	F6	Indole-3-propionic Acid	28821	Aromatic amino acid derivative
1	F7	L-Tryptophan	29600	Aromatic amino acid derivative
1	F8	Indole-3-pyruvic Acid	29876	Aromatic amino acid derivative
1	F9	4-Hydroxyhippuric Acid	30115	Aromatic amino acid derivative
1	F10	<i>Unused</i>		
1	F11	Indole-3-carboxylic Acid	31166	Aromatic amino acid derivative
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 8

Plate	Well	Contents	Item Number	Class
1	G1	<i>Unused</i>		
1	G2	Histamine	33828	Aromatic amino acid derivative
1	G3	Indole-3-acetamide	34374	Aromatic amino acid derivative
1	G4	N-(3-Phenylpropionyl)glycine	34392	Aromatic amino acid derivative
1	G5	Salicylic Acid	34813	Aromatic amino acid derivative
1	G6	<i>Unused</i>		
1	G7	<i>Unused</i>		
1	G8	<i>Unused</i>		
1	G9	<i>Unused</i>		
1	G10	<i>Unused</i>		
1	G11	<i>Unused</i>		
1	G12	<i>Unused</i>		
1	H1	<i>Unused</i>		
1	H2	<i>Unused</i>		
1	H3	<i>Unused</i>		
1	H4	<i>Unused</i>		
1	H5	<i>Unused</i>		
1	H6	<i>Unused</i>		
1	H7	<i>Unused</i>		
1	H8	<i>Unused</i>		
1	H9	<i>Unused</i>		
1	H10	<i>Unused</i>		
1	H11	<i>Unused</i>		
1	H12	<i>Unused</i>		
2	A1	<i>Unused</i>		
2	A2	(±)-Enterolactone	10112	Aromatic
2	A3	Ellagic Acid	10569	Aromatic
2	A4	Delphinidin (chloride)	11012	Aromatic
2	A5	Coumestrol	11730	Aromatic
2	A6	(-)-Epicatechin	11807	Aromatic
2	A7	Gallic Acid	11846	Aromatic
2	A8	Kaempferol	11852	Aromatic
2	A9	Safrole	13894	Aromatic
2	A10	Glycitein	14162	Aromatic
2	A11	Naringenin	14173	Aromatic
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 8

Plate	Well	Contents	Item Number	Class
2	B1	<i>Unused</i>		
2	B2	N-Acetylserotonin	14535	Aromatic
2	B3	Cyanidin (chloride)	14803	Aromatic
2	B4	Protocatechuic Acid	14916	Aromatic
2	B5	(±)-Salsolinol (hydrochloride)	15597	Aromatic
2	B6	Isorhamnetin	16496	Aromatic
2	B7	1-Methylnicotinamide (chloride)	16604	Aromatic
2	B8	Naringin	17923	Aromatic
2	B9	4-Aminobenzoic Acid (sodium salt)	18659	Aromatic
2	B10	Dihydroresveratrol	19651	Aromatic
2	B11	Malvidin (chloride)	19752	Aromatic
2	B12	<i>Unused</i>		
2	C1	<i>Unused</i>		
2	C2	Pelargonidin (chloride)	19753	Aromatic
2	C3	Peonidin (chloride)	19754	Aromatic
2	C4	Petunidin (chloride)	19755	Aromatic
2	C5	Ferulic Acid	19871	Aromatic
2	C6	Pyrogallol	20347	Aromatic
2	C7	(±)-Equol 4'-sulfate (sodium salt)	21440	Aromatic
2	C8	N-Methylnicotinamide	21850	Aromatic
2	C9	Urolithin A	22607	Aromatic
2	C10	(+)-Gallocatechin	25125	Aromatic
2	C11	Sinapic Acid	25457	Aromatic
2	C12	<i>Unused</i>		
2	D1	<i>Unused</i>		
2	D2	2S-Eriodictyol	26803	Aromatic
2	D3	Homovanillic Acid	27307	Aromatic
2	D4	4-Hydroxybenzoic Acid	27468	Aromatic
2	D5	Gentisic Acid	27615	Aromatic
2	D6	(±)-Catechin (hydrate)	29175	Aromatic
2	D7	<i>p</i> -Cresol sulfate (potassium salt)	29504	Aromatic
2	D8	Salicylic Acid	32991	Aromatic
2	D9	4-Aminosalicylic Acid	33449	Aromatic
2	D10	Urolithin B	33757	Aromatic
2	D11	Prunin	34034	Aromatic
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



Plate	Well	Contents	Item Number	Class
2	E1	<i>Unused</i>		
2	E2	Nicotinic Acid	34096	Aromatic
2	E3	Catechol	34791	Aromatic
2	E4	5-Aminosalicylic Acid	70265	Aromatic
2	E5	Guaiacol	70430	Aromatic
2	E6	Caffeic Acid	70602	Aromatic
2	E7	<i>trans</i> -Resveratrol	70675	Aromatic
2	E8	Luteolin	10004161	Aromatic
2	E9	Daidzein	10005166	Aromatic
2	E10	Genistein	10005167	Aromatic
2	E11	Quercetin (hydrate)	10005169	Aromatic
2	E12	<i>Unused</i>		
2	F1	<i>Unused</i>		
2	F2	(S)-Equol	10010173	Aromatic
2	F3	Apigenin	10010275	Aromatic
2	F4	Myricetin	10012600	Aromatic
2	F5	<i>Unused</i>		
2	F6	<i>Unused</i>		
2	F7	<i>Unused</i>		
2	F8	<i>Unused</i>		
2	F9	Piperine	11750	Dietary alkaloid
2	F10	<i>Unused</i>		
2	F11	Quinine	23958	Dietary alkaloid
2	F12	<i>Unused</i>		
2	G1	<i>Unused</i>		
2	G2	Sulforaphane	10496	Dietary terpene, lignan, and glucosinolate
2	G3	Sesamol	11754	Dietary terpene, lignan, and glucosinolate
2	G4	β -Carotene	16837	Dietary terpene, lignan, and glucosinolate
2	G5	β -Caryophyllene	21572	Dietary terpene, lignan, and glucosinolate
2	G6	β -Myrcene	21720	Dietary terpene, lignan, and glucosinolate
2	G7	(\pm)-Limonene	25773	Dietary terpene, lignan, and glucosinolate
2	G8	(+)-Pinoresinol	27700	Dietary terpene, lignan, and glucosinolate
2	G9	Sesamin	70310	Dietary terpene, lignan, and glucosinolate
2	G10	Matairesinol	10005174	Dietary terpene, lignan, and glucosinolate
2	G11	Glucoraphanin (potassium salt)	10009445	Dietary terpene, lignan, and glucosinolate
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



Plate	Well	Contents	Item Number	Class
2	H1	Unused		
2	H2	Unused		
2	H3	Azelaic Acid	23977	Dioic acid
2	H4	3-Methylglutaric Acid	34357	Dioic acid
2	H5	CMPF	10007133	Dioic acid
2	H6	Unused		
2	H7	Unused		
2	H8	Unused		
2	H9	Unused		
2	H10	Unused		
2	H11	Unused		
2	H12	Unused		
3	A1	Unused		
3	A2	(R)-3-hydroxy Myristic Acid	16871	Lipid
3	A3	Commendamide	18421	Lipid
3	A4	3-hydroxy Palmitic Acid	19934	Lipid
3	A5	1-Oleoyl-2-hydroxy-sn-glycero-3-PG (sodium salt)	20960	Lipid
3	A6	1-Palmitoyl-2-hydroxy-sn-glycero-3-PG (sodium salt)	20961	Lipid
3	A7	2-hydroxy Palmitic Acid	22679	Lipid
3	A8	3-hydroxy Stearic Acid	22692	Lipid
3	A9	2-hydroxy Lauric Acid	24586	Lipid
3	A10	2-hydroxy Stearic Acid	24590	Lipid
3	A11	3-hydroxy Lauric Acid	24642	Lipid
3	A12	Unused		
3	B1	Unused		
3	B2	11-methyl Lauric Acid methyl ester	24809	Lipid
3	B3	12-methyl Tridecanoic Acid methyl ester	24810	Lipid
3	B4	13-methyl Myristic Acid	24811	Lipid
3	B5	14-methyl Pentadecanoic Acid methyl ester	24813	Lipid
3	B6	15-methyl Palmitic Acid	24814	Lipid
3	B7	12-methyl Myristic Acid	24817	Lipid
3	B8	14-methyl Palmitic Acid	24820	Lipid
3	B9	Unused		
3	B10	17-methyl Stearic Acid	26766	Lipid
3	B11	3-hydroxy Heptadecanoic Acid	28735	Lipid
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 8

Plate	Well	Contents	Item Number	Class
3	C1	Unused		
3	C2	2-hydroxy Myristic Acid	90390	Lipid
3	C3	Palmitoleic Acid	10009871	Lipid
3	C4	Unused		
3	C5	Unused		
3	C6	Unused		
3	C7	1,3,7-Trimethyluric Acid	16949	Nucleotide metabolite
3	C8	1,7-Dimethyluric Acid	19584	Nucleotide metabolite
3	C9	Allantoin	30204	Nucleotide metabolite
3	C10	3-Methylxanthine	34288	Nucleotide metabolite
3	C11	7-Methylxanthine	34348	Nucleotide metabolite
3	C12	Unused		
3	D1	Unused		
3	D2	N-Formyl-Met-Leu-Phe	21495	Peptide
3	D3	Muramyl Dipeptide	30866	Peptidoglycan component
3	D4	N-Acetylmuramic Acid	31330	Peptidoglycan component
3	D5	Unused		
3	D6	Unused		
3	D7	Unused		
3	D8	Unused		
3	D9	Unused		
3	D10	Unused		
3	D11	Unused		
3	D12	Unused		
3	E1	Unused		
3	E2	Unused		
3	E3	(±)-β-Hydroxybutyrate*	14148	SCFA/Small aliphatic acid
3	E4	Unused		
3	E5	Octanoic Acid*	33674	SCFA/Small aliphatic acid
3	E6	α-Ketoglutaric Acid*	33693	SCFA/Small aliphatic acid
3	E7	β-Hydroxyisovaleric Acid*	34030	SCFA/Small aliphatic acid
3	E8	α-Ketoisocaproic Acid*	34749	SCFA/Small aliphatic acid
3	E9	Unused		
3	E10	Unused		
3	E11	Unused		
3	E12	Unused		

*The item formulation for this item is: A 1 M solution in DMSO.

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 8

Plate	Well	Contents	Item Number	Class
3	F1	Unused		
3	F2	C18 Ceramide (d18:1/18:0)	19556	Sphingolipid
3	F3	Unused		
3	F4	Unused		
3	F5	Unused		
3	F6	C17 Ceramide (d18:1/17:0)	22532	Sphingolipid
3	F7	Unused		
3	F8	Unused		
3	F9	Unused		
3	F10	Sphingosine (d17:1)	10007902	Sphingolipid
3	F11	Sphingosine (d18:1)	10007907	Sphingolipid
3	F12	Unused		
3	G1	Unused		
3	G2	Sphinganine (d18:0)	10007945	Sphingolipid
3	G3	Unused		
3	G4	Unused		
3	G5	Unused		
3	G6	Unused		
3	G7	Unused		
3	G8	Unused		
3	G9	Unused		
3	G10	Unused		
3	G11	Coprostanol	26764	Sterol
3	G12	Unused		
3	H1	Unused		
3	H2	Unused		
3	H3	Unused		
3	H4	Unused		
3	H5	Unused		
3	H6	Unused		
3	H7	Unused		
3	H8	Unused		
3	H9	Unused		
3	H10	Unused		
3	H11	Unused		
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