

PRODUCT INFORMATION



Synthetic Cannabinoid Analytical Standards Panel 3

Item No. 38770

Synonym: Synthetic CB Analytical Standards Panel 3
Supplied As: 100 µg per well, each component at a nominal 1 mg/ml concentration
Storage: -20°C
Stability: ≥4 years

Laboratory Procedures

Each tube contains one synthetic cannabinoid supplied at a nominal 1 mg/ml concentration with 100 µl provided. Before opening, warm each plate to room temperature while upright to allow the solution within each tube to settle to the bottom. Use proper Personal Protective Equipment (PPE) when handling the tube wells. The solution can be used as-is, but if a lower concentration and higher volume is preferred, an example dilution procedure follows:

1. Uncap each tube to be used. Add 400 µl of the solvent corresponding to the diluent column on the table below to each tube. This will create 500 µl of a 200 µg/ml solution of each analyte. *NOTE: Tube capacity is 700 µl. No more than 500 µl of solvent should be added.*
2. Recap the tube(s) and place the plate on a plate mixer or vortex mixer. Plates and/or tube wells should be well mixed or vortexed to ensure a homogenous solution within the tube.
3. Carefully remove the cap(s) for analysis.

Description

Synthetic Cannabinoid Analytical Standards Panel 3 contains 204 analytical reference standards categorized as synthetic cannabinoids and synthetic cannabinoid metabolites. This curated collection encompasses the wave of synthetic cannabinoid compounds that became relevant to the forensic community from 2017 through 2024. For the generation of synthetic cannabinoids and synthetic cannabinoid metabolites preceding this timespan, see Synthetic Cannabinoid Analytical Standards Panel 1 (Item No. 38768) or Synthetic Cannabinoid Analytical Standards Panel 2 (Item No. 38769). The compounds in Synthetic Cannabinoid Analytical Standards Panel 3 are supplied at a nominal 1 mg/ml concentration with 100 µl provided in each tube of a 96-well tube rack for rapid screening or library building.

Please review the product table, below, for a full list of targets. The Synthetic Cannabinoid Analytical Standards Panel 3 contains compounds regulated as Schedule I compounds in the United States and is regulated as a Schedule I item. This product is intended for research and forensic applications.

Panel Contents

Plate	Well	Contents	Item Number	Diluent	Molecular Formula	CAS Number
1	A1	<i>Unused</i>				
1	A2	CUMYL-PINACA	17207	Acetonitrile	C ₂₂ H ₂₇ N ₃ O	1400742-15-5
1	A3	ADB-FUBICA	18756	Acetonitrile	C ₂₂ H ₂₄ FN ₃ O ₂	2748161-90-0
1	A4	MEP-FUBICA	21653	Acetonitrile	C ₂₂ H ₂₃ FN ₂ O ₃	2749985-70-2
1	A5	ADB-5'Br-PINACA 3,3-dimethylbutanoic acid metabolite	42007	Acetonitrile	C ₁₉ H ₂₆ BrN ₃ O ₃	N/A
1	A6	5-fluoro ADB metabolite 7	20178	Acetonitrile	C ₁₉ H ₂₆ FN ₃ O ₃	2365471-39-0
1	A7	ADB-5'F-BUTINACA	37761	Acetonitrile	C ₁₈ H ₂₅ FN ₄ O ₂	N/A
1	A8	5-fluoro MDMB-PICA metabolite 7	22095	Acetonitrile	C ₂₀ H ₂₇ FN ₂ O ₃	2569223-22-7
1	A9	MEP-CHMICA	21652	Acetonitrile	C ₂₂ H ₃₀ N ₂ O ₃	2749985-65-5
1	A10	NAPIE	22632	Acetonitrile	C ₂₅ H ₂₅ NO	2748289-69-0
1	A11	NAMIE	22631	Acetonitrile	C ₂₁ H ₁₇ NO	1638677-49-2
1	A12	<i>Unused</i>				

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, 10/02/2025

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	Diluent	Molecular Formula	CAS Number
1	B1	<i>Unused</i>				
1	B2	7'-methoxy NABUTIE	22630	Acetonitrile	C ₂₅ H ₂₅ NO ₂	1438278-55-7
1	B3	5-fluoro CUMYL-PINACA N-pentanoic acid metabolite	31286	Acetonitrile	C ₂₂ H ₂₅ N ₃ O ₃	N/A
1	B4	4-fluoro MDMB-BUTINACA 3-carboxyindazole metabolite	27819	Acetonitrile	C ₁₂ H ₁₃ FN ₂ O ₂	2027536-90-7
1	B5	MDMB-INACA 3,3-dimethylbutanoic acid metabolite	38151	Acetonitrile	C ₁₄ H ₁₇ N ₃ O ₃	1189552-82-6
1	B6	5-fluoro EDMB-PINACA	23005	Acetonitrile	C ₂₁ H ₃₀ FN ₃ O ₃	2504100-69-8
1	B7	5-fluoro- <i>tert</i> -Butylbenzyl-PINACA	23351	Acetonitrile	C ₂₄ H ₃₀ FN ₃ O	2749299-04-3
1	B8	(S)-5-fluoro ADB	23631	Acetonitrile	C ₂₀ H ₂₈ FN ₃ O ₃	1971007-89-2
1	B9	Ethylphenethyl-FUBICA	23764	Acetonitrile	C ₂₆ H ₂₅ FN ₂ O	2749394-77-0
1	B10	EADB-FUBINACA	23345	Acetonitrile	C ₂₃ H ₂₇ FN ₄ O ₂	2749394-81-6
1	B11	CUMYL-PeGACLONE	22696	Acetonitrile	C ₂₅ H ₂₈ N ₂ O	2160555-55-3
1	B12	<i>Unused</i>				
1	C1	<i>Unused</i>				
1	C2	AZEFUBIM	23653	Acetonitrile	C ₁₉ H ₁₇ FN ₂ O	2748592-03-0
1	C3	MEP-FUBINACA	21654	Acetonitrile	C ₂₁ H ₂₂ FN ₃ O ₃	2749985-46-2
1	C4	Ethylbenzyl-CYBINACA	24431	Acetonitrile	C ₂₂ H ₂₄ N ₄ O	2749618-88-8
1	C5	4-fluoro MDMB-BUTICA butanoic acid metabolite	31106	Acetonitrile	C ₁₉ H ₂₅ FN ₂ O ₃	N/A
1	C6	MDMB-PINACA	39730	Acetonitrile	C ₂₀ H ₂₉ N ₃ O ₃	1971007-99-4
1	C7	Methyl 1-(4-fluorobenzyl)-1H-indazole-3-carboxylate	24855	Acetonitrile	C ₁₆ H ₁₃ FN ₂ O ₂	1185287-50-6
1	C8	CH-PIATA N-pentanoic acid metabolite	37603	Acetonitrile	C ₂₁ H ₂₈ N ₂ O ₃	N/A
1	C9	5-fluoro 7-APAICA	23021	Acetonitrile	C ₂₃ H ₃₀ FN ₃ O	2682867-58-7
1	C10	ACHMINACA	25006	Acetonitrile	C ₂₅ H ₃₃ N ₃ O	1400742-33-7
1	C11	5-fluoro 7-QUPAIC	23011	Acetonitrile	C ₂₂ H ₂₀ FN ₃ O ₂	2748300-92-5
1	C12	<i>Unused</i>				
1	D1	<i>Unused</i>				
1	D2	4-fluoro ADB	42153	Acetonitrile	C ₂₀ H ₂₈ FN ₃ O ₃	2365471-11-8
1	D3	5-fluoro ADB metabolite 2	20173	Acetonitrile	C ₂₀ H ₂₉ N ₃ O ₄	2642091-65-2
1	D4	FUBIMINA N-(5-hydroxypentyl) metabolite	24468	Acetonitrile	C ₂₃ H ₂₂ N ₂ O ₂	2748343-95-3
1	D5	MDMB-5'Br-BUTINACA	37698	Acetonitrile	C ₁₉ H ₂₆ BrN ₃ O ₃	N/A
1	D6	MMB-FUBICA metabolite 3	24619	Acetonitrile	C ₂₁ H ₂₁ FN ₂ O ₃	2693397-44-1
1	D7	5-bromo APINACA	25209	Acetonitrile	C ₂₃ H ₃₀ BrN ₃ O	2160555-51-9
1	D8	MMB022	25906	Acetonitrile	C ₂₀ H ₂₆ N ₂ O ₃	2659308-31-1
1	D9	Isobutyl 1-pentyl-1H-indazole-3-carboxylate	25880	Acetonitrile	C ₁₇ H ₂₄ N ₂ O ₂	2748624-88-4
1	D10	MDMB-4en-PINACA	26097	Acetonitrile	C ₂₀ H ₂₇ N ₃ O ₃	2504100-70-1
1	D11	5-fluoro phenyl-PICA	23924	Acetonitrile	C ₂₀ H ₂₁ FN ₂ O	1776086-01-1
1	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	Diluent	Molecular Formula	CAS Number
1	E1	<i>Unused</i>				
1	E2	4-chloro CUMYL-PINACA	26132	Acetonitrile	C ₂₂ H ₂₆ ClN ₃ O	1631074-65-1
1	E3	4-fluoro-CUMYL-5-fluoro-PICA	26131	Acetonitrile	C ₂₃ H ₂₆ F ₂ N ₂ O	1631074-52-6
1	E4	5-fluoro MPP-PICA	25916	Acetonitrile	C ₂₄ H ₂₇ FN ₂ O ₃	2682867-54-3
1	E5	4-fluoro-CUMYL-5-fluoro-PINACA	26134	Acetonitrile	C ₂₂ H ₂₅ F ₂ N ₃ O	1631074-53-7
1	E6	5-fluoro MDMA-7-PAICA	25286	Acetonitrile	C ₂₀ H ₂₈ FN ₃ O ₃	2377403-81-9
1	E7	4-fluoro MDMA-BUTINACA	26645	Acetonitrile	C ₁₉ H ₂₆ FN ₃ O ₃	2390036-46-9
1	E8	APP-BUTINACA	26905	Acetonitrile	C ₂₁ H ₂₄ N ₄ O ₂	2682867-56-5
1	E9	MDMA-INACA	27248	Acetonitrile	C ₁₅ H ₁₉ N ₃ O ₃	2709672-58-0
1	E10	4-fluoro MDMA-BUTINACA butanoic acid metabolite	26856	Acetonitrile	C ₁₈ H ₂₄ FN ₃ O ₃	N/A
1	E11	MDMA-5Me-INACA	39658	Acetonitrile	C ₁₆ H ₂₁ N ₃ O ₃	N/A
1	E12	<i>Unused</i>				
1	F1	<i>Unused</i>				
1	F2	4-fluoro MDMA-BUTINACA N-(4-hydroxybutyl) metabolite	26857	Acetonitrile	C ₁₉ H ₂₇ N ₃ O ₄	N/A
1	F3	4-fluoro MDMA-BUTINACA 2'-indazole isomer	27442	Acetonitrile	C ₁₉ H ₂₆ FN ₃ O ₃	2706356-20-7
1	F4	AB-PINACA 3-carboxyindazole metabolite	27732	Acetonitrile	C ₁₃ H ₁₆ N ₂ O ₂	1283576-17-9
1	F5	5-fluoro AB-PINACA 3-carboxyindazole metabolite	27733	Acetonitrile	C ₁₃ H ₁₅ FN ₂ O ₂	1535166-43-8
1	F6	CUMYL-NB-MeGACLONE	32793	Acetonitrile	C ₂₈ H ₃₀ N ₂ O	N/A
1	F7	AB-MDMSBA	41967	Acetonitrile	C ₁₅ H ₂₃ N ₃ O ₄ S	N/A
1	F8	6-fluoro MDA 19	40803	Acetonitrile	C ₂₁ H ₂₂ FN ₃ O ₂	N/A
1	F9	MDMA-CHMICA metabolite M2	19316	Acetonitrile	C ₂₂ H ₃₀ N ₂ O ₃	2460730-12-3
1	F10	5-fluoro MDMA-PICA metabolite 8	27586	Acetonitrile	C ₁₉ H ₂₄ N ₂ O ₅	N/A
1	F11	4-cyano MDMA-BUTINACA	26507	Acetonitrile	C ₂₀ H ₂₆ N ₄ O ₃	1185888-30-5
1	F12	<i>Unused</i>				
1	G1	<i>Unused</i>				
1	G2	2-fluoro QMPSB	27325	Acetonitrile	C ₂₂ H ₂₀ F ₂ N ₂ O ₄ S	2707165-48-6
1	G3	3TMS-CUMYL-PRINACA	40509	Acetonitrile	C ₂₃ H ₃₁ N ₃ OSi	N/A
1	G4	5-fluoro MDMA-PICA metabolite 2	22393	Acetonitrile	C ₂₁ H ₃₀ N ₂ O ₄	N/A
1	G5	MDMA-3en-BUTINACA	28508	Acetonitrile	C ₁₉ H ₂₅ N ₃ O ₃	N/A
1	G6	MDMA-4en-PICA	28541	Acetonitrile	C ₂₁ H ₂₈ N ₂ O ₃	2659308-32-2
1	G7	MDMA-4en-PINACA butanoic acid metabolite	28740	Acetonitrile	C ₁₉ H ₂₅ N ₃ O ₃	N/A
1	G8	2-fluoro AMB	37319	Acetonitrile	C ₁₉ H ₂₆ FN ₃ O ₃	N/A
1	G9	5-fluoro MDMA-PICA metabolite 9	27587	Acetonitrile	C ₂₁ H ₂₉ FN ₂ O ₄	N/A
1	G10	MDMA-BUTINACA	28739	Acetonitrile	C ₁₉ H ₂₇ N ₃ O ₃	3039541-81-3
1	G11	3TMS-ADB-PRINACA	38844	Acetonitrile	C ₂₀ H ₃₂ N ₄ O ₂ Si	N/A
1	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	Diluent	Molecular Formula	CAS Number
1	H1	<i>Unused</i>				
1	H2	MMB-ICA	39951	Acetonitrile	C ₁₅ H ₁₈ N ₂ O ₃	1188516-52-0
1	H3	5-fluoro ADB 2'-indazole isomer	28962	Acetonitrile	C ₂₀ H ₂₈ FN ₃ O ₃	2712429-78-0
1	H4	MDMB-MINACA	39659	Acetonitrile	C ₁₆ H ₂₁ N ₃ O ₃	N/A
1	H5	PTI-3	28238	Acetonitrile	C ₂₁ H ₂₈ FN ₃ OS	N/A
1	H6	NE-CHMIMO	29301	Acetonitrile	C ₂₆ H ₂₅ NO	1373876-11-9
1	H7	ADB-BUTINACA	29350	Acetonitrile	C ₁₈ H ₂₆ N ₄ O ₂	2682867-55-4
1	H8	5-bromo AMB-PICA	29752	Acetonitrile	C ₂₀ H ₂₇ BrN ₂ O ₃	N/A
1	H9	MMB-4en-PINACA	29750	Acetonitrile	C ₁₉ H ₂₅ N ₃ O ₃	2659308-41-3
1	H10	MMB-4en-PINACA butanoic acid metabolite	30301	Acetonitrile	C ₁₈ H ₂₃ N ₃ O ₃	N/A
1	H11	AB-5Br-INACA	39953	Acetonitrile	C ₁₃ H ₁₅ BrN ₄ O ₂	N/A
1	H12	<i>Unused</i>				
2	A1	<i>Unused</i>				
2	A2	MDMB-ICA	27249	Acetonitrile	C ₁₆ H ₂₀ N ₂ O ₃	2741575-87-9
2	A3	5-chloro MDMB-PICA	23438	Acetonitrile	C ₂₁ H ₂₉ ClN ₂ O ₃	N/A
2	A4	5-chloro AMB-PICA	29751	Acetonitrile	C ₂₀ H ₂₇ ClN ₂ O ₃	N/A
2	A5	BUTINACA	40096	Acetonitrile	C ₁₂ H ₁₅ N ₃ O	3036872-65-5
2	A6	Methyl 2-(1H-indazole-3-carboxamido)-3-methylbutanoate	25580	Acetonitrile	C ₁₄ H ₁₇ N ₃ O ₃	1563540-10-2
2	A7	4-fluoro ABUTINACA	30724	Acetonitrile	C ₂₂ H ₂₈ FN ₃ O	1445580-39-1
2	A8	5-fluoro EMB-PICA	30769	Acetonitrile	C ₂₁ H ₂₉ FN ₂ O ₃	2648861-83-8
2	A9	ADB-BUTINACA N-butanoic acid metabolite	30294	Acetonitrile	C ₁₈ H ₂₄ N ₄ O ₄	N/A
2	A10	MDMB-BUTICA	40205	Acetonitrile	C ₂₀ H ₂₈ N ₂ O ₃	N/A
2	A11	5-fluoro EDMB-PICA	30725	Acetonitrile	C ₂₂ H ₃₁ FN ₂ O ₃	2666934-54-7
2	A12	<i>Unused</i>				
2	B1	<i>Unused</i>				
2	B2	MDMB-4en-PICA butanoic acid metabolite	28943	Acetonitrile	C ₂₀ H ₂₆ N ₂ O ₃	N/A
2	B3	MDMB-3en-BUTINACA butanoic acid metabolite	28942	Acetonitrile	C ₁₈ H ₂₃ N ₃ O ₃	N/A
2	B4	4-fluoro MDMB-BUTICA	31075	Acetonitrile	C ₂₀ H ₂₇ FN ₂ O ₃	2682867-53-2
2	B5	CUMYL-PIPETINACA (hydrochloride)	27331	Acetonitrile	C ₂₄ H ₃₀ N ₄ O • HCl	N/A
2	B6	iPDMB-FUBINACA	28930	Acetonitrile	C ₂₄ H ₂₈ FN ₃ O ₃	N/A
2	B7	AB-INACA	39400	Acetonitrile	C ₁₃ H ₁₆ N ₄ O ₂	1887742-41-7
2	B8	5-fluoro MMB-P7AICA	31189	Acetonitrile	C ₁₉ H ₂₆ FN ₃ O ₃	2755646-64-9
2	B9	4-cyano CUMYL-BUT7AICA	22818	Acetonitrile	C ₂₂ H ₂₄ N ₄ O	2160555-53-1
2	B10	MPP-PICA	31268	Acetonitrile	C ₂₄ H ₂₈ N ₂ O ₃	N/A
2	B11	Benzyl-4-cyano BUTINACA	31313	Acetonitrile	C ₂₀ H ₂₀ N ₄ O	N/A
2	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



Plate	Well	Contents	Item Number	Diluent	Molecular Formula	CAS Number
2	C1	<i>Unused</i>				
2	C2	CUMYL-CBMINACA	31325	Acetonitrile	C ₂₂ H ₂₅ N ₃ O	N/A
2	C3	MDMB-PICA	39729	Acetonitrile	C ₂₁ H ₃₀ N ₂ O ₃	1971007-98-3
2	C4	MMB-CHM7AICA	31188	Acetonitrile	C ₂₁ H ₂₉ N ₃ O ₃	2749438-69-3
2	C5	4-fluoro ABUTINACA N-butanoic acid metabolite	30749	Acetonitrile	C ₂₂ H ₂₇ N ₃ O ₃	N/A
2	C6	4-cyano ADB-BUTINACA	26506	Acetonitrile	C ₁₉ H ₂₅ N ₅ O ₂	N/A
2	C7	AMB butanoic acid metabolite	30302	Acetonitrile	C ₁₈ H ₂₅ N ₃ O ₃	1890250-16-4
2	C8	5-fluoro MMB-PICA butanoic acid metabolite	31267	Acetonitrile	C ₁₉ H ₂₅ FN ₂ O ₃	2693397-43-0
2	C9	4-fluoro MDMB-BUTICA N-(4-hydroxybutyl) metabolite	31541	Acetonitrile	C ₂₀ H ₂₈ N ₂ O ₄	N/A
2	C10	4-cyano CUMYL-BUTCZCA	24429	Acetonitrile	C ₂₇ H ₂₇ N ₃ O	2740821-84-3
2	C11	MDMB-CHM7AICA	31186	Acetonitrile	C ₂₂ H ₃₁ N ₃ O ₃	2741576-83-8
2	C12	<i>Unused</i>				
2	D1	<i>Unused</i>				
2	D2	5-fluoro EMB-PICA N-(5-hydroxypentyl) metabolite	31540	Acetonitrile	C ₂₁ H ₃₀ N ₂ O ₄	N/A
2	D3	MDMB-FUB7AICA	31187	Acetonitrile	C ₂₂ H ₂₄ FN ₃ O ₃	2751267-51-1
2	D4	4-fluoro EMB-BUTICA	31298	Acetonitrile	C ₂₀ H ₂₇ FN ₂ O ₃	N/A
2	D5	THQ-PINAC	27081	Acetonitrile	C ₂₂ H ₂₅ N ₃ O	N/A
2	D6	ADB-BUTINACA N-(4-hydroxybutyl) metabolite	30295	Acetonitrile	C ₁₈ H ₂₆ N ₄ O ₃	N/A
2	D7	4-chloro MDMB-BUTICA	31805	Acetonitrile	C ₂₀ H ₂₇ ClN ₂ O ₃	N/A
2	D8	APP-BUTINACA phenylpropanoic acid metabolite	31279	Acetonitrile	C ₂₁ H ₂₃ N ₃ O ₃	N/A
2	D9	4-chloro MDMB-BUTINACA	42183	Acetonitrile	C ₁₉ H ₂₆ ClN ₃ O ₃	N/A
2	D10	MMB-4en-PICA butanoic acid metabolite	31275	Acetonitrile	C ₁₉ H ₂₄ N ₂ O ₃	N/A
2	D11	CUMYL-CB-MeGACLONE	31427	Acetonitrile	C ₂₅ H ₂₆ N ₂ O	2806439-13-2
2	D12	<i>Unused</i>				
2	E1	<i>Unused</i>				
2	E2	4-fluoro MDMB-BUTINACA N-(butanoic acid) 3,3-dimethylbutanoic acid metabolite	31643	Acetonitrile	C ₁₈ H ₂₃ N ₃ O ₅	N/A
2	E3	AFUB7AICA	31185	Acetonitrile	C ₂₅ H ₂₆ FN ₃ O	2810271-54-4
2	E4	ADB-BUTINAATA	36014	Acetonitrile	C ₁₉ H ₂₈ N ₄ O ₂	N/A
2	E5	ADB-4en-PINACA	33205	Acetonitrile	C ₁₉ H ₂₆ N ₄ O ₂	2666932-44-9
2	E6	4-fluoro EDMB-BUTINACA	33374	Acetonitrile	C ₂₀ H ₂₈ FN ₃ O ₃	N/A
2	E7	JWH 184	32794	Acetonitrile	C ₂₅ H ₂₇ N	619294-37-0
2	E8	5-fluoro CUMYL-PeGACLONE N-pentanoic acid metabolite	31284	Acetonitrile	C ₂₅ H ₂₆ N ₂ O ₃	2835243-96-2
2	E9	5-fluoro CUMYL-PeGACLONE N-(5-hydroxypentyl) metabolite	31283	Acetonitrile	C ₂₅ H ₂₈ N ₂ O ₂	2399452-31-2
2	E10	4-cyano MMB-BUTINACA	33334	Acetonitrile	C ₁₉ H ₂₄ N ₄ O ₃	2659308-48-0
2	E11	4-fluoro ABUTINACA N-(4-hydroxybutyl) metabolite	30750	Acetonitrile	C ₂₂ H ₂₉ N ₃ O ₂	N/A
2	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



Plate	Well	Contents	Item Number	Diluent	Molecular Formula	CAS Number
2	F1	<i>Unused</i>				
2	F2	MMB-FUBGACLONE	30770	Acetonitrile	C ₂₄ H ₂₃ FN ₂ O ₃	2797662-15-6
2	F3	4-cyano MMB-BUTINACA butanoic acid metabolite	33668	Acetonitrile	C ₁₈ H ₂₂ N ₄ O ₃	N/A
2	F4	ADB-PHETINACA	33194	Acetonitrile	C ₂₂ H ₂₆ N ₄ O ₂	N/A
2	F5	MDMB-5'Br-4en-PINACA	39488	Acetonitrile	C ₂₀ H ₂₆ BrN ₃ O ₃	N/A
2	F6	AMP-4en-PINACA	33818	Acetonitrile	C ₁₇ H ₂₂ N ₄ O ₂	N/A
2	F7	ADB-HEXINACA	33820	Acetonitrile	C ₂₀ H ₃₀ N ₄ O ₂	N/A
2	F8	EDMB-PINACA	33663	Acetonitrile	C ₂₁ H ₃₁ N ₃ O ₃	2666934-55-8
2	F9	NMDMSB	38886	Acetonitrile	C ₂₀ H ₁₉ NO ₄ S	N/A
2	F10	CHO-4'Me-5'Br-FUBOXPYRA	37468	Acetonitrile	C ₂₀ H ₂₂ BrFN ₂ O ₂	N/A
2	F11	4-cyano MMB-BUTINACA N-butanoic acid metabolite	33835	Acetonitrile	C ₁₈ H ₂₃ N ₃ O ₅	N/A
2	F12	<i>Unused</i>				
2	G1	<i>Unused</i>				
2	G2	BZO-POXIZID	34795	Acetonitrile	C ₂₀ H ₂₁ N ₃ O ₂	1048973-64-3
2	G3	MDMB-BENZICA	35135	Acetonitrile	C ₂₃ H ₂₆ N ₂ O ₃	N/A
2	G4	5-fluoro BZO-POXIZID	34586	Acetonitrile	C ₂₀ H ₂₀ FN ₃ O ₂	N/A
2	G5	ADB-P7AICA	34352	Acetonitrile	C ₁₉ H ₂₈ N ₄ O ₂	2366273-07-4
2	G6	4-cyano MMB-BUTINACA N-butanoic acid 3-methylbutanoic acid metabolite	33836	Acetonitrile	C ₁₇ H ₂₁ N ₃ O ₅	N/A
2	G7	CUMYL-NBMINACA	34908	Acetonitrile	C ₂₅ H ₂₉ N ₃ O	1631074-60-6
2	G8	AEP-CHMINACA	23706	Acetonitrile	C ₂₀ H ₂₈ N ₄ O ₂	N/A
2	G9	AEP-PINACA	23707	Acetonitrile	C ₁₈ H ₂₆ N ₄ O ₂	N/A
2	G10	4-fluoro AB-BUTINACA	37657	Acetonitrile	C ₁₇ H ₂₃ FN ₄ O ₂	1864889-46-2
2	G11	ADB-FUBIATA	35059	Acetonitrile	C ₂₃ H ₂₆ FN ₃ O ₂	2938025-73-9
2	G12	<i>Unused</i>				
2	H1	<i>Unused</i>				
2	H2	BZO-4en-POXIZID	35503	Acetonitrile	C ₂₀ H ₁₉ N ₃ O ₂	N/A
2	H3	2-(1-(4-Fluorobenzyl)-1H-indol-3-yl)acetic Acid	36015	Acetonitrile	C ₁₇ H ₁₄ FNO ₂	176204-51-6
2	H4	MDA-19 N-(5-hydroxyhexyl) metabolite	36039	Acetonitrile	C ₂₁ H ₂₃ N ₃ O ₃	N/A
2	H5	ADB-FUBIATA 3,3-dimethylbutanoic acid metabolite	35222	Acetonitrile	C ₂₃ H ₂₅ FN ₂ O ₃	N/A
2	H6	JWH 181	38120	Acetonitrile	C ₂₈ H ₃₁ NO	824960-03-4
2	H7	ADB-HEXINACA N-(6-hydroxyhexyl) metabolite	35566	Acetonitrile	C ₂₀ H ₃₀ N ₄ O ₃	N/A
2	H8	ADB-HEXINACA N-hexanoic acid metabolite	35773	Acetonitrile	C ₂₀ H ₂₈ N ₄ O ₄	N/A
2	H9	5-fluoro ADB-P7AICA	34353	Acetonitrile	C ₁₉ H ₂₇ FN ₄ O ₂	N/A
2	H10	MDA 19	10563	Acetonitrile	C ₂₁ H ₂₃ N ₃ O ₂	1048973-47-2
2	H11	3,5-ADB-4en-PFUPPYCA	36141	Acetonitrile	C ₂₁ H ₂₇ FN ₄ O ₂	N/A
2	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



Plate	Well	Contents	Item Number	Diluent	Molecular Formula	CAS Number
3	A1	<i>Unused</i>				
3	A2	EDMB-4en-PINACA	36444	Acetonitrile	C ₂₁ H ₂₉ N ₃ O ₃	N/A
3	A3	MDMB-5Br-INACA	36757	Acetonitrile	C ₁₅ H ₁₈ BrN ₃ O ₃	N/A
3	A4	5,3-ADB-4en-PFUPPYCA	36146	Acetonitrile	C ₂₁ H ₂₇ FN ₄ O ₂	N/A
3	A5	ADB-HEXINACA 3,3-dimethylbutanoic acid metabolite	35567	Acetonitrile	C ₂₀ H ₂₉ N ₃ O ₃	N/A
3	A6	5-fluoro ADB-5'Br-PINACA	37963	Acetonitrile	C ₁₉ H ₂₆ BrFN ₄ O ₂	N/A
3	A7	CH-PIATA	36558	Acetonitrile	C ₂₁ H ₃₀ N ₂ O	N/A
3	A8	ADB-5'Br-BUTINACA 3,3-dimethyl butanoic acid metabolite	37607	Acetonitrile	C ₁₈ H ₂₄ BrN ₃ O ₃	N/A
3	A9	MPP-BUTINACA	36852	Acetonitrile	C ₂₂ H ₂₅ N ₃ O ₃	N/A
3	A10	5-fluoro AMB metabolite 6	37058	Acetonitrile	C ₁₈ H ₂₃ N ₃ O ₅	1890250-15-3
3	A11	MMB-5Br-INACA	36883	Acetonitrile	C ₁₄ H ₁₆ BrN ₃ O ₃	N/A
3	A12	<i>Unused</i>				
3	B1	<i>Unused</i>				
3	B2	AFUBIATA	36992	Acetonitrile	C ₂₇ H ₂₉ FN ₂ O	2985275-14-5
3	B3	CUMYL-CHSINACA	37057	Acetonitrile	C ₂₃ H ₂₇ N ₃ O ₃ S	N/A
3	B4	ADB-FUBINAATA	36853	Acetonitrile	C ₂₂ H ₂₅ FN ₄ O ₂	N/A
3	B5	BZO-HEPOXIZID	36946	Acetonitrile	C ₂₂ H ₂₅ N ₃ O ₂	N/A
3	B6	<i>Unused</i>				
3	B7	<i>Unused</i>				
3	B8	<i>Unused</i>				
3	B9	<i>Unused</i>				
3	B10	<i>Unused</i>				
3	B11	<i>Unused</i>				
3	B12	<i>Unused</i>				
3	C1	<i>Unused</i>				
3	C2	<i>Unused</i>				
3	C3	<i>Unused</i>				
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

PRODUCT INFORMATION



Plate	Well	Contents	Item Number	Diluent	Molecular Formula	CAS Number
3	D1	Unused				
3	D2	Unused				
3	D3	Unused				
3	D4	Unused				
3	D5	Unused				
3	D6	Unused				
2	D7	Unused				
2	D8	Unused				
2	D9	Unused				
3	D10	Unused				
3	D11	Unused				
3	D12	Unused				
3	E1	Unused				
3	E2	Unused				
3	E3	Unused				
3	E4	Unused				
3	E5	Unused				
3	E6	Unused				
3	E7	Unused				
3	E8	Unused				
3	E9	Unused				
3	E10	Unused				
3	E11	Unused				
3	E12	Unused				
3	F1	Unused				
3	F2	CUMYL-PICA N-pentanoic acid metabolite	20232	Methanol	C ₂₃ H ₂₆ N ₂ O ₃	2748623-20-1
3	F3	FUBIMINA N-pentanoic acid metabolite	24469	Methanol	C ₂₃ H ₂₀ N ₂ O ₃	2748624-01-1
3	F4	4-fluoro MDMB-BUTINACA 3-carboxy-2'-indazole metabolite	27820	Methanol	C ₁₂ H ₁₃ FN ₂ O ₂	2709837-98-7
3	F5	5-fluoro CUMYL-PeGACLONE	25435	Methanol	C ₂₅ H ₂₇ FN ₂ O	2377403-49-9
3	F6	4-cyano CUMYL-BUTINACA metabolite 10	26399	Methanol	C ₂₁ H ₂₃ N ₃ O ₃	N/A
3	F7	4-fluoro MDMB-BUTINACA 2'-isomer butanoic acid metabolite	28964	Methanol	C ₁₈ H ₂₄ FN ₃ O ₃	2713210-84-3
3	F8	5-fluoro ADB 2'-indazole isomer butanoic acid metabolite	28963	Methanol	C ₁₉ H ₂₆ FN ₃ O ₃	2711867-70-6
3	F9	2-(2-Chlorophenyl)-1-(1H-indol-3-yl)ethanone	39498	Methanol	C ₁₆ H ₁₂ ClNO	1096914-22-5
3	F10	CUMYL-CBMICA	30207	Methanol	C ₂₃ H ₂₆ N ₂ O	2571070-88-5
3	F11	MDMB-BUTINACA butanoic acid metabolite	28939	Methanol	C ₁₈ H ₂₅ N ₃ O ₃	3039541-82-4
3	F12	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



Plate	Well	Contents	Item Number	Diluent	Molecular Formula	CAS Number
3	G1	<i>Unused</i>				
3	G2	5-fluoro AB-7-PAICA	26646	Methanol	C ₁₈ H ₂₅ FN ₄ O ₂	2682867-57-6
3	G3	4-fluoro MDMA-BUTINACA N-butanoic acid metabolite	28941	Methanol	C ₁₉ H ₂₅ N ₃ O ₅	N/A
3	G4	Ethylbenzyl-PeGACLONE	9003353	Methanol	C ₂₅ H ₂₈ N ₂ O	2805900-54-1
2	G5	BZO-CHMOXIZID	34956	Methanol	C ₂₂ H ₂₃ N ₃ O ₂	1048973-67-6
2	G6	AFUB7AICA 7'-azaindole isomer	32989	Methanol	C ₂₅ H ₂₆ FN ₃ O	2809146-29-8
2	G7	ADB-5Br-INACA	36263	Methanol	C ₁₄ H ₁₇ BrN ₄ O ₂	N/A
3	G8	ADB-5'Br-BUTINACA	36770	Methanol	C ₁₈ H ₂₅ BrN ₄ O ₂	N/A
3	G9	CH-FUBIATA	37054	Methanol	C ₂₃ H ₂₅ FN ₂ O	922038-77-5
3	G10	5-fluoro MPP-PICA phenylpropionic acid metabolite	31271	Methanol	C ₂₃ H ₂₅ FN ₂ O ₃	N/A
3	G11	CUMYL-INACA	37055	Methanol	C ₁₇ H ₁₇ N ₃ O	1631075-21-2
3	G12	<i>Unused</i>				
3	H1	<i>Unused</i>				
3	H2	ADB-5'Br-DECINACA	37721	Methanol	C ₂₄ H ₃₇ BrN ₄ O ₂	N/A
3	H3	ADB-5'Br-PINACA	37758	Methanol	C ₁₉ H ₂₇ BrN ₄ O ₂	N/A
3	H4	ADB-5'Br-4en-PINACA	37741	Methanol	C ₁₉ H ₂₅ BrN ₄ O ₂	N/A
3	H5	ADB-INACA	37729	Methanol	C ₁₄ H ₁₈ N ₄ O ₂	1887742-42-8
3	H6	ADB-IATA	37053	Methanol	C ₁₆ H ₂₁ N ₃ O ₂	N/A
3	H7	CH-HEXIATA	38283	Methanol	C ₂₂ H ₃₂ N ₂ O	N/A
3	H8	CHM-FUBIATA	38205	Methanol	C ₂₄ H ₂₇ FN ₂ O	922092-52-2
3	H9	5-fluoro MDMA-7-PAICA butanoic acid metabolite	28693	Methanol	C ₁₉ H ₂₆ FN ₃ O ₃	2712863-53-9
3	H10	4-cyano AB-BUTICA	35192	Methanol	C ₁₉ H ₂₄ N ₄ O ₂	N/A
3	H11	MN-18 N-(5-hydroxypentyl) metabolite	35278	Methanol	C ₂₃ H ₂₃ N ₃ O ₂	2748591-74-2
3	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