

CERTIFICATE of ANALYSIS



Synthetic Cannabinoid Analytical Standards Panel 3

For R&D purposes only. Not intended for human or animal use.

Item No.:	38770
Batch No.:	0808202
Size:	100 µg/well
Expiry Date:	19SEP2029 (valid from date of certification)
Storage & Handling:	Store upright and unopened at -20°C. Warm to room temperature prior to opening.
Safety:	Danger. Review the Safety Data Sheets before use.

Description

This product consists of three 96-well Matrix plates containing individual synthetic cannabinoids, metabolites, and precursors. The individual well locations for each unique standard are listed in the table below.

Traceability Statement

This panel was prepared from individual standards that have been verified through standard quality testing and are traceable via individual item and batch numbers. Individual certificates of analysis are available upon request. All compounds used to make this panel were used within their valid expiry dates. Stability testing of this panel is ongoing to ensure continuous quality and traceability of the materials.

Quality Testing

A single panel was pulled randomly from the batch and mass analysis was performed to verify the contents of each vial position.

Test	Specification	Meets Specification
Analyte Detected, MS	± 0.5 amu	Yes

Individual results are reported in the table on pages 2-13.

Approval:

Title: Manager, Analytical Chemistry

Certification Date: 19SEP2025

Cayman Chemical certifies that this product meets the specifications stated in this certificate and warrants this product to meet the stated acceptance criteria through the expiry date when stored unopened as recommended.

CERTIFICATE of ANALYSIS



Kit Contents Plate 1

Well	Item Name	Molecular Formula (free analyte)	Expected [M+H] ⁺ m/z	Observed [M+H] ⁺ m/z
A1	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
A2	CUMYL-PINACA	C ₂₂ H ₂₇ N ₃ O	350.2	350.2
A3	ADB-FUBICA	C ₂₂ H ₂₄ FN ₃ O ₂	382.2	382.3
A4	MEP-FUBICA	C ₂₂ H ₂₃ FN ₂ O ₃	383.2	383.1
A5	ADB-5'Br-PINACA 3,3-dimethylbutanoic acid metabolite	C ₁₉ H ₂₆ BrN ₃ O ₃	424.1	424.1
A6	5-fluoro ADB metabolite 7	C ₁₉ H ₂₆ FN ₃ O ₃	364.2	363.9
A7	ADB-5'F-BUTINACA	C ₁₈ H ₂₅ FN ₄ O ₂	349.2	349.1
A8	5-fluoro MDMA-PICA metabolite 7	C ₂₀ H ₂₇ FN ₂ O ₃	363.2	363.1
A9	MEP-CHMICA	C ₂₂ H ₃₀ N ₂ O ₃	371.2	371.1
A10	NAPIE	C ₂₅ H ₂₅ NO	356.2	356.1
A11	NAMIE	C ₂₁ H ₁₇ NO	300.1	300.1
A12	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
B1	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
B2	7'-methoxy NABUTIE	C ₂₅ H ₂₅ NO ₂	372.2	372
B3	5-fluoro CUMYL-PINACA N-pentanoic acid metabolite	C ₂₂ H ₂₅ N ₃ O ₃	380.2	380.1
B4	4-fluoro MDMA-BUTINACA 3-carboxyindazole metabolite	C ₁₂ H ₁₃ FN ₂ O ₂	237.1	237.2
B5	MDMA-INACA 3,3-dimethylbutanoic acid metabolite	C ₁₄ H ₁₇ N ₃ O ₃	276.1	275.9
B6	5-fluoro MDMA-PINACA	C ₂₁ H ₃₀ FN ₃ O ₃	392.2	392
B7	5-fluoro- <i>tert</i> -Butylbenzyl-PINACA	C ₂₄ H ₃₀ FN ₃ O	396.2	396
B8	(S)-5-fluoro ADB	C ₂₀ H ₂₈ FN ₃ O ₃	378.2	378.2
B9	Ethylphenethyl-FUBICA	C ₂₆ H ₂₅ FN ₂ O	401.2	401.1
B10	EADB-FUBINACA	C ₂₃ H ₂₇ FN ₄ O ₂	411.2	411.2
B11	CUMYL-PeGACLONE	C ₂₅ H ₂₈ N ₂ O	373.2	373.1
B12	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<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

techserv@caymanchem.com
www.caymanchem.com

CERTIFICATE of ANALYSIS



Kit Contents Plate 1 Continued

Well	Item Name	Molecular Formula (free analyte)	Expected [M+H] ⁺ m/z	Observed [M+H] ⁺ m/z
C1	Unused	Unused	Unused	Unused
C2	AZEFUBIM	C ₁₉ H ₁₇ FN ₂ O	309.1	309.1
C3	MEP-FUBINACA	C ₂₁ H ₂₂ FN ₃ O ₃	384.2	384.4
C4	Ethylbenzyl-CYBINACA	C ₂₂ H ₂₄ N ₄ O	361.2	361.2
C5	4-fluoro MDMA-BUTICA butanoic acid metabolite	C ₁₉ H ₂₅ FN ₂ O ₃	349.2	349.1
C6	MDMA-PINACA	C ₂₀ H ₂₉ N ₃ O ₃	360.2	360.1
C7	Methyl 1-(4-fluorobenzyl)-1H-indazole-3-carboxylate	C ₁₆ H ₁₃ FN ₂ O ₂	285.1	285.2
C8	CH-PIATA N-pentanoic acid metabolite	C ₂₁ H ₂₈ N ₂ O ₃	357.2	357.2
C9	5-fluoro 7-APAICA	C ₂₃ H ₃₀ FN ₃ O	384.2	384.1
C10	ACHMINACA	C ₂₅ H ₃₃ N ₃ O	392.3	392.2
C11	5-fluoro 7-QUPAIC	C ₂₂ H ₂₀ FN ₃ O ₂	378.2	377.9
C12	Unused	Unused	Unused	Unused
D1	Unused	Unused	Unused	Unused
D2	4-fluoro ADB	C ₂₀ H ₂₈ FN ₃ O ₃	378.2	378
D3	5-fluoro ADB metabolite 2	C ₂₀ H ₂₉ N ₃ O ₄	376.2	376.1
D4	FUBIMINA N-(5-hydroxypentyl) metabolite	C ₂₃ H ₂₂ N ₂ O ₂	359.2	358.8
D5	MDMA-5'Br-BUTINACA	C ₁₉ H ₂₆ BrN ₃ O ₃	424.1	424.1
D6	MMB-FUBICA metabolite 3	C ₂₁ H ₂₁ FN ₂ O ₃	369.2	369.1
D7	5-bromo APINACA	C ₂₃ H ₃₀ BrN ₃ O	444.2	444.1
D8	MMB022	C ₂₀ H ₂₆ N ₂ O ₃	343.2	343.1
D9	Isobutyl 1-pentyl-1H-indazole-3-carboxylate	C ₁₇ H ₂₄ N ₂ O ₂	289.2	289.2
D10	MDMA-4en-PINACA	C ₂₀ H ₂₇ N ₃ O ₃	358.2	358.1
D11	5-fluoro phenyl-PICA	C ₂₀ H ₂₁ FN ₂ O	325.2	325.1
D12	Unused	Unused	Unused	Unused

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640

techserv@caymanchem.com
www.caymanchem.com

CERTIFICATE of ANALYSIS



Kit Contents Plate 1 Continued

Well	Item Name	Molecular Formula (free analyte)	Expected [M+H] ⁺ m/z	Observed [M+H] ⁺ m/z
E1	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
E2	4-chloro CUMYL-PINACA	C ₂₂ H ₂₆ ClN ₃ O	384.2	383.9
E3	4-fluoro-CUMYL-5-fluoro-PICA	C ₂₃ H ₂₆ F ₂ N ₂ O	385.2	385.2
E4	5-fluoro MPP-PICA	C ₂₄ H ₂₇ FN ₂ O ₃	411.2	411.2
E5	4-fluoro-CUMYL-5-fluoro-PINACA	C ₂₂ H ₂₅ F ₂ N ₃ O	386.2	386.1
E6	5-fluoro MDMA-7-PAICA	C ₂₀ H ₂₈ FN ₃ O ₃	378.2	378.1
E7	4-fluoro MDMA-BUTINACA	C ₁₉ H ₂₆ FN ₃ O ₃	364.2	364.1
E8	APP-BUTINACA	C ₂₁ H ₂₄ N ₄ O ₂	365.2	365.2
E9	MDMA-INACA	C ₁₅ H ₁₉ N ₃ O ₃	290.1	290.2
E10	4-fluoro MDMA-BUTINACA butanoic acid metabolite	C ₁₈ H ₂₄ FN ₃ O ₃	350.2	350.1
E11	MDMA-5Me-INACA	C ₁₆ H ₂₁ N ₃ O ₃	304.2	303.9
E12	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
F1	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
F2	4-fluoro MDMA-BUTINACA N-(4-hydroxybutyl) metabolite	C ₁₉ H ₂₇ N ₃ O ₄	362.2	362
F3	4-fluoro MDMA-BUTINACA 2'-indazole isomer	C ₁₉ H ₂₆ FN ₃ O ₃	364.2	364.1
F4	AB-PINACA 3-carboxyindazole metabolite	C ₁₃ H ₁₆ N ₂ O ₂	233.1	233.2
F5	5-fluoro AB-PINACA 3-carboxyindazole metabolite	C ₁₃ H ₁₅ FN ₂ O ₂	251.1	251.2
F6	CUMYL-NB-MeGACLONE	C ₂₈ H ₃₀ N ₂ O	411.2	411
F7	AB-MDMSBA	C ₁₅ H ₂₃ N ₃ O ₄ S	342.1	342.1
F8	6-fluoro MDA 19	C ₂₁ H ₂₂ FN ₃ O ₂	368.2	368.1
F9	MDMA-CHMICA metabolite M2	C ₂₂ H ₃₀ N ₂ O ₃	371.2	371.2
F10	5-fluoro MDMA-PICA metabolite 8	C ₁₉ H ₂₄ N ₂ O ₅	361.2	361.1
F11	4-cyano MDMA-BUTINACA	C ₂₀ H ₂₆ N ₄ O ₃	371.2	371.1
F12	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<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

techserv@caymanchem.com
www.caymanchem.com

CERTIFICATE of ANALYSIS



Kit Contents Plate 1 Continued

Well	Item Name	Molecular Formula (free analyte)	Expected [M+H] ⁺ m/z	Observed [M+H] ⁺ m/z
G1	Unused	Unused	Unused	Unused
G2	2-fluoro QMPSB	C ₂₂ H ₂₀ F ₂ N ₂ O ₄ S	447.1	447.1
G3	3TMS-CUMYL-PRINACA	C ₂₃ H ₃₁ N ₃ OSi	394.2	394.1
G4	5-fluoro MDMA-PICA metabolite 2	C ₂₁ H ₃₀ N ₂ O ₄	375.2	375.3
G5	MDMA-3en-BUTINACA	C ₁₉ H ₂₅ N ₃ O ₃	344.2	344.1
G6	MDMA-4en-PICA	C ₂₁ H ₂₈ N ₂ O ₃	357.2	357.2
G7	MDMA-4en-PINACA butanoic acid metabolite	C ₁₉ H ₂₅ N ₃ O ₃	344.2	344.1
G8	2-fluoro AMB	C ₁₉ H ₂₆ FN ₃ O ₃	364.2	364.1
G9	5-fluoro MDMA-PICA metabolite 9	C ₂₁ H ₂₉ FN ₂ O ₄	393.2	393.2
G10	MDMA-BUTINACA	C ₁₉ H ₂₇ N ₃ O ₃	346.2	346.2
G11	3TMS-ADB-PRINACA	C ₂₀ H ₃₂ N ₄ O ₂ Si	389.2	389.1
G12	Unused	Unused	Unused	Unused
H1	Unused	Unused	Unused	Unused
H2	MMB-ICA	C ₁₅ H ₁₈ N ₂ O ₃	275.1	275.2
H3	5-fluoro ADB 2'-indazole isomer	C ₂₀ H ₂₈ FN ₃ O ₃	378.2	378.1
H4	MDMA-MINACA	C ₁₆ H ₂₁ N ₃ O ₃	304.2	304.2
H5	PTI-3	C ₂₁ H ₂₈ FN ₃ OS	390.2	390.2
H6	NE-CHMIMO	C ₂₆ H ₂₅ NO	368.2	368
H7	ADB-BUTINACA	C ₁₈ H ₂₆ N ₄ O ₂	331.2	331.1
H8	5-bromo AMB-PICA	C ₂₀ H ₂₇ BrN ₂ O ₃	423.1	423.2
H9	MMB-4en-PINACA	C ₁₉ H ₂₅ N ₃ O ₃	344.2	344.1
H10	MMB-4en-PINACA butanoic acid metabolite	C ₁₈ H ₂₃ N ₃ O ₃	330.2	330.1
H11	AB-5Br-INACA	C ₁₃ H ₁₅ BrN ₄ O ₂	339	339.2
H12	Unused	Unused	Unused	Unused

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640

techserv@caymanchem.com
www.caymanchem.com

CERTIFICATE of ANALYSIS



Kit Contents Plate 2

Well	Item Name	Molecular Formula (free analyte)	Expected [M+H] ⁺ m/z	Observed [M+H] ⁺ m/z
A1	Unused	Unused	Unused	Unused
A2	MDMB-ICA	C ₁₆ H ₂₀ N ₂ O ₃	289.2	289.2
A3	5-chloro MDMB-PICA	C ₂₁ H ₂₉ ClN ₂ O ₃	393.2	392.9
A4	5-chloro AMB-PICA	C ₂₀ H ₂₇ ClN ₂ O ₃	379.2	379.1
A5	BUTINACA	C ₁₂ H ₁₅ N ₃ O	218.1	218.2
A6	Methyl 2-(1H-indazole-3-carboxamido)-3-methylbutanoate	C ₁₄ H ₁₇ N ₃ O ₃	276.1	276.1
A7	4-fluoro ABUTINACA	C ₂₂ H ₂₈ FN ₃ O	370.2	370.1
A8	5-fluoro EMB-PICA	C ₂₁ H ₂₉ FN ₂ O ₃	377.2	377.1
A9	ADB-BUTINACA N-butanoic acid metabolite	C ₁₈ H ₂₄ N ₄ O ₄	361.2	361.1
A10	MDMB-BUTICA	C ₂₀ H ₂₈ N ₂ O ₃	345.2	345.2
A11	5-fluoro EDMB-PICA	C ₂₂ H ₃₁ FN ₂ O ₃	391.2	391.2
A12	Unused	Unused	Unused	Unused
B1	Unused	Unused	Unused	Unused
B2	MDMB-4en-PICA butanoic acid metabolite	C ₂₀ H ₂₆ N ₂ O ₃	343.2	343.1
B3	MDMB-3en-BUTINACA butanoic acid metabolite	C ₁₈ H ₂₃ N ₃ O ₃	330.2	330.1
B4	4-fluoro MDMB-BUTICA	C ₂₀ H ₂₇ FN ₂ O ₃	363.2	363.3
B5	CUMYL-PIPETINACA (hydrochloride)	C ₂₄ H ₃₀ N ₄ O	391.2	391.2
B6	iPDMB-FUBINACA	C ₂₄ H ₂₈ FN ₃ O ₃	426.2	426.1
B7	AB-INACA	C ₁₃ H ₁₆ N ₄ O ₂	261.1	261.2
B8	5-fluoro MMB-P7AICA	C ₁₉ H ₂₆ FN ₃ O ₃	364.2	364
B9	4-cyano CUMYL-BUT7AICA	C ₂₂ H ₂₄ N ₄ O	361.2	361.1
B10	MPP-PICA	C ₂₄ H ₂₈ N ₂ O ₃	393.2	393.2
B11	Benzyl-4-cyano BUTINACA	C ₂₀ H ₂₀ N ₄ O	333.2	333.1
B12	Unused	Unused	Unused	Unused

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640

techserv@caymanchem.com
www.caymanchem.com

CERTIFICATE of ANALYSIS



Kit Contents Plate 2 Continued

Well	Item Name	Molecular Formula (free analyte)	Expected [M+H] ⁺ m/z	Observed [M+H] ⁺ m/z
C1	Unused	Unused	Unused	Unused
C2	CUMYL-CBMINACA	C ₂₂ H ₂₅ N ₃ O	348.2	348.1
C3	MDMB-PICA	C ₂₁ H ₃₀ N ₂ O ₃	359.2	359.2
C4	MMB-CHM7AICA	C ₂₁ H ₂₉ N ₃ O ₃	372.2	372.1
C5	4-fluoro ABUTINACA N-butanoic acid metabolite	C ₂₂ H ₂₇ N ₃ O ₃	382.2	382.2
C6	4-cyano ADB-BUTINACA	C ₁₉ H ₂₅ N ₅ O ₂	356.2	356
C7	AMB butanoic acid metabolite	C ₁₈ H ₂₅ N ₃ O ₃	332.2	331.8
C8	5-fluoro MMB-PICA butanoic acid metabolite	C ₁₉ H ₂₅ FN ₂ O ₃	349.2	349.2
C9	4-fluoro MDMB-BUTICA N-(4-hydroxybutyl) metabolite	C ₂₀ H ₂₈ N ₂ O ₄	361.2	361.1
C10	4-cyano CUMYL-BUTCZCA	C ₂₇ H ₂₇ N ₃ O	410.2	410.2
C11	MDMB-CHM7AICA	C ₂₂ H ₃₁ N ₃ O ₃	386.2	386.2
C12	Unused	Unused	Unused	Unused
D1	Unused	Unused	Unused	Unused
D2	5-fluoro EMB-PICA N-(5-hydroxypentyl) metabolite	C ₂₁ H ₃₀ N ₂ O ₄	375.2	375.2
D3	MDMB-FUB7AICA	C ₂₂ H ₂₄ FN ₃ O ₃	398.2	398.2
D4	4-fluoro EMB-BUTICA	C ₂₀ H ₂₇ FN ₂ O ₃	363.2	363.1
D5	THQ-PINACA	C ₂₂ H ₂₅ N ₃ O	348.2	347.9
D6	ADB-BUTINACA N-(4-hydroxybutyl) metabolite	C ₁₈ H ₂₆ N ₄ O ₃	347.2	347.1
D7	4-chloro MDMB-BUTICA	C ₂₀ H ₂₇ ClN ₂ O ₃	379.2	379
D8	APP-BUTINACA phenylpropanoic acid metabolite	C ₂₁ H ₂₃ N ₃ O ₃	366.2	366.1
D9	4-chloro MDMB-BUTINACA	C ₁₉ H ₂₆ ClN ₃ O ₃	380.2	380.1
D10	MMB-4en-PICA butanoic acid metabolite	C ₁₉ H ₂₄ N ₂ O ₃	329.2	329
D11	CUMYL-CB-MeGACLONE	C ₂₅ H ₂₆ N ₂ O	371.2	371.2
D12	Unused	Unused	Unused	Unused

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640

techserv@caymanchem.com
www.caymanchem.com

CERTIFICATE of ANALYSIS



Kit Contents Plate 2 Continued

Well	Item Name	Molecular Formula (free analyte)	Expected [M+H] ⁺ m/z	Observed [M+H] ⁺ m/z
E1	Unused	Unused	Unused	Unused
E2	4-fluoro MDMB-BUTINACA N-(butanoic acid) 3,3-dimethylbutanoic acid metabolite	C ₁₈ H ₂₃ N ₃ O ₅	362.2	362
E3	AFUB7AICA	C ₂₅ H ₂₆ FN ₃ O	404.2	404.1
E4	ADB-BUTINAATA	C ₁₉ H ₂₈ N ₄ O ₂	345.2	345.2
E5	ADB-4en-PINACA	C ₁₉ H ₂₆ N ₄ O ₂	343.2	343.2
E6	4-fluoro EDMB-BUTINACA	C ₂₀ H ₂₈ FN ₃ O ₃	378.2	378.2
E7	JWH 184	C ₂₅ H ₂₇ N	342.2	342.2
E8	5-fluoro CUMYL-PeGACLONE N-pentanoic acid metabolite	C ₂₅ H ₂₆ N ₂ O ₃	403.2	403.2
E9	5-fluoro CUMYL-PeGACLONE N-(5-hydroxypentyl) metabolite	C ₂₅ H ₂₈ N ₂ O ₂	389.2	389.1
E10	4-cyano MMB-BUTINACA	C ₁₉ H ₂₄ N ₄ O ₃	357.2	357.1
E11	4-fluoro ABUTINACA N-(4-hydroxybutyl) metabolite	C ₂₂ H ₂₉ N ₃ O ₂	368.2	368
E12	Unused	Unused	Unused	Unused
F1	Unused	Unused	Unused	Unused
F2	MMB-FUBGACLONE	C ₂₄ H ₂₃ FN ₂ O ₃	407.2	407.1
F3	4-cyano MMB-BUTINACA butanoic acid metabolite	C ₁₈ H ₂₂ N ₄ O ₃	343.2	343
F4	ADB-PHETINACA	C ₂₂ H ₂₆ N ₄ O ₂	379.2	379.2
F5	MDMB-5'Br-4en-PINACA	C ₂₀ H ₂₆ BrN ₃ O ₃	436.1	436
F6	AMP-4en-PINACA	C ₁₇ H ₂₂ N ₄ O ₂	315.2	315.1
F7	ADB-HEXINACA	C ₂₀ H ₃₀ N ₄ O ₂	359.2	359.2
F8	EDMB-PINACA	C ₂₁ H ₃₁ N ₃ O ₃	374.2	374.2
F9	NMDMSB	C ₂₀ H ₁₉ NO ₄ S	370.1	370.1
F10	CHO-4'Me-5'Br-FUBOXPYRA	C ₂₀ H ₂₂ BrFN ₂ O ₂	421.1	421
F11	4-cyano MMB-BUTINACA N-butanoic acid metabolite	C ₁₈ H ₂₃ N ₃ O ₅	362.2	362.1
F12	Unused	Unused	Unused	Unused

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640

techserv@caymanchem.com
www.caymanchem.com

CERTIFICATE of ANALYSIS



Kit Contents Plate 2 Continued

Well	Item Name	Molecular Formula (free analyte)	Expected [M+H] ⁺ m/z	Observed [M+H] ⁺ m/z
G1	Unused	Unused	Unused	Unused
G2	BZO-POXIZID	C ₂₀ H ₂₁ N ₃ O ₂	336.2	336.1
G3	MDMB-BENZICA	C ₂₃ H ₂₆ N ₂ O ₃	379.2	379.2
G4	5-fluoro BZO-POXIZID	C ₂₀ H ₂₀ FN ₃ O ₂	354.2	354
G5	ADB-P7AICA	C ₁₉ H ₂₈ N ₄ O ₂	345.2	345.2
G6	4-cyano MMB-BUTINACA N-butanolic acid 3-methylbutanoic acid metabolite	C ₁₇ H ₂₁ N ₃ O ₅	348.2	348.1
G7	CUMYL-NBMINACA	C ₂₅ H ₂₉ N ₃ O	388.2	388.2
G8	AEP-CHMINACA	C ₂₀ H ₂₈ N ₄ O ₂	357.2	357.1
G9	AEP-PINACA	C ₁₈ H ₂₆ N ₄ O ₂	331.2	331.2
G10	4-fluoro AB-BUTINACA	C ₁₇ H ₂₃ FN ₄ O ₂	335.2	335.1
G11	ADB-FUBIATA	C ₂₃ H ₂₆ FN ₃ O ₂	396.2	396.1
G12	Unused	Unused	Unused	Unused
H1	Unused	Unused	Unused	Unused
H2	BZO-4en-POXIZID	C ₂₀ H ₁₉ N ₃ O ₂	334.2	334.1
H3	2-(1-(4-Fluorobenzyl)-1H-indol-3-yl)acetic Acid	C ₁₇ H ₁₄ FNO ₂	284.1	284.1
H4	MDA-19 N-(5-hydroxyhexyl) metabolite	C ₂₁ H ₂₃ N ₃ O ₃	366.2	366.2
H5	ADB-FUBIATA 3,3-dimethylbutanoic acid metabolite	C ₂₃ H ₂₅ FN ₃ O ₃	397.2	397.2
H6	JWH 181	C ₂₈ H ₃₁ NO	398.2	398.2
H7	ADB-HEXINACA N-(6-hydroxyhexyl) metabolite	C ₂₀ H ₃₀ N ₄ O ₃	375.2	375.2
H8	ADB-HEXINACA N-hexanoic acid metabolite	C ₂₀ H ₂₈ N ₄ O ₄	389.2	389.2
H9	5-fluoro ADB-P7AICA	C ₁₉ H ₂₇ FN ₄ O ₂	363.2	363.2
H10	MDA 19	C ₂₁ H ₂₃ N ₃ O ₂	350.2	350.1
H11	3,5-ADB-4en-PFUPPYCA	C ₂₁ H ₂₇ FN ₄ O ₂	387.2	387
H12	Unused	Unused	Unused	Unused

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640

techserv@caymanchem.com
www.caymanchem.com

CERTIFICATE of ANALYSIS



Kit Contents Plate 3

Well	Item Name	Molecular Formula (free analyte)	Expected [M+H] ⁺ m/z	Observed [M+H] ⁺ m/z
A1	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
A2	EDMB-4en-PINACA	C ₂₁ H ₂₉ N ₃ O ₃	372.2	372.2
A3	MDMB-5Br-INACA	C ₁₅ H ₁₈ BrN ₃ O ₃	368.1	368
A4	5,3-ADB-4en-PFUPPYCA	C ₂₁ H ₂₇ FN ₄ O ₂	387.2	387.2
A5	ADB-HEXINACA 3,3-dimethylbutanoic acid metabolite	C ₂₀ H ₂₉ N ₃ O ₃	360.2	360.1
A6	5-fluoro ADB-5'Br-PINACA	C ₁₉ H ₂₆ BrFN ₄ O ₂	441.1	441.1
A7	CH-PIATA	C ₂₁ H ₃₀ N ₂ O	327.2	327
A8	ADB-5'Br-BUTINACA 3,3-dimethyl butanoic acid metabolite	C ₁₈ H ₂₄ BrN ₃ O ₃	410.1	410
A9	MPP-BUTINACA	C ₂₂ H ₂₅ N ₃ O ₃	380.2	380
A10	5-fluoro AMB metabolite 6	C ₁₈ H ₂₃ N ₃ O ₅	362.2	362
A11	MMB-5Br-INACA	C ₁₄ H ₁₆ BrN ₃ O ₃	354	353.9
A12	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
B1	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
B2	AFUBIATA	C ₂₇ H ₂₉ FN ₂ O	417.2	417
B3	CUMYL-CHSINACA	C ₂₃ H ₂₇ N ₃ O ₃ S	426.2	426.1
B4	ADB-FUBINAATA	C ₂₂ H ₂₅ FN ₄ O ₂	397.2	397.1
B5	BZO-HEPOXIZID	C ₂₂ H ₂₅ N ₃ O ₂	364.2	364.2
B6	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
B7	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
B8	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
B9	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
B10	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
B11	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
B12	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<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

techserv@caymanchem.com
www.caymanchem.com

CERTIFICATE of ANALYSIS



Kit Contents Plate 3 Continued

Well	Item Name	Molecular Formula (free analyte)	Expected [M+H] ⁺ m/z	Observed [M+H] ⁺ m/z
C1	Unused	Unused	Unused	Unused
C2	Unused	Unused	Unused	Unused
C3	Unused	Unused	Unused	Unused
C4	Unused	Unused	Unused	Unused
C5	Unused	Unused	Unused	Unused
C6	Unused	Unused	Unused	Unused
C7	Unused	Unused	Unused	Unused
C8	Unused	Unused	Unused	Unused
C9	Unused	Unused	Unused	Unused
C10	Unused	Unused	Unused	Unused
C11	Unused	Unused	Unused	Unused
C12	Unused	Unused	Unused	Unused
D1	Unused	Unused	Unused	Unused
D2	Unused	Unused	Unused	Unused
D3	Unused	Unused	Unused	Unused
D4	Unused	Unused	Unused	Unused
D5	Unused	Unused	Unused	Unused
D6	Unused	Unused	Unused	Unused
D7	Unused	Unused	Unused	Unused
D8	Unused	Unused	Unused	Unused
D9	Unused	Unused	Unused	Unused
D10	Unused	Unused	Unused	Unused
D11	Unused	Unused	Unused	Unused
D12	Unused	Unused	Unused	Unused

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640

techserv@caymanchem.com
www.caymanchem.com

CERTIFICATE of ANALYSIS



Kit Contents Plate 3 Continued

Well	Item Name	Molecular Formula (free analyte)	Expected [M+H] ⁺ m/z	Observed [M+H] ⁺ m/z
E1	Unused	Unused	Unused	Unused
E2	Unused	Unused	Unused	Unused
E3	Unused	Unused	Unused	Unused
E4	Unused	Unused	Unused	Unused
E5	Unused	Unused	Unused	Unused
E6	Unused	Unused	Unused	Unused
E7	Unused	Unused	Unused	Unused
E8	Unused	Unused	Unused	Unused
E9	Unused	Unused	Unused	Unused
E10	Unused	Unused	Unused	Unused
E11	Unused	Unused	Unused	Unused
E12	Unused	Unused	Unused	Unused
F1	Unused	Unused	Unused	Unused
F2	CUMYL-PICA N-pentanoic acid metabolite	C ₂₃ H ₂₆ N ₂ O ₃	379.2	379.2
F3	FUBIMINA N-pentanoic acid metabolite	C ₂₃ H ₂₀ N ₂ O ₃	373.2	373
F4	4-fluoro MDMA-BUTINACA 3-carboxy-2'-indazole metabolite	C ₁₂ H ₁₃ FN ₂ O ₂	237.1	237.1
F5	5-fluoro CUMYL-PeGACLONE	C ₂₅ H ₂₇ FN ₂ O	391.2	391.1
F6	4-cyano CUMYL-BUTINACA metabolite 10	C ₂₁ H ₂₃ N ₃ O ₃	366.2	366
F7	4-fluoro MDMA-BUTINACA 2'-isomer butanoic acid metabolite	C ₁₈ H ₂₄ FN ₃ O ₃	350.2	350.1
F8	5-fluoro ADB 2'-indazole isomer butanoic acid metabolite	C ₁₉ H ₂₆ FN ₃ O ₃	364.2	364.1
F9	2-(2-Chlorophenyl)-1-(1H-indol-3-yl)ethanone	C ₁₆ H ₁₂ ClNO	270.1	270.1
F10	CUMYL-CBMICA	C ₂₃ H ₂₆ N ₂ O	347.2	347.1
F11	MDMA-BUTINACA butanoic acid metabolite	C ₁₈ H ₂₅ N ₃ O ₃	332.2	332.1
F12	Unused	Unused	Unused	Unused

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640

techserv@caymanchem.com
www.caymanchem.com

CERTIFICATE of ANALYSIS



Kit Contents Plate 3 Continued

Well	Item Name	Molecular Formula (free analyte)	Expected [M+H] ⁺ m/z	Observed [M+H] ⁺ m/z
G1	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
G2	5-fluoro AB-7-PAICA	C ₁₈ H ₂₅ FN ₄ O ₂	349.2	349.1
G3	4-fluoro MDMA-BUTINACA N-butanoic acid metabolite	C ₁₉ H ₂₅ N ₃ O ₅	376.2	376
G4	Ethylbenzyl-PeGACLONE	C ₂₅ H ₂₈ N ₂ O	373.2	373.2
G5	BZO-CHMOXIZID	C ₂₂ H ₂₃ N ₃ O ₂	362.2	362
G6	AFUB7AICA 7'-azaindole isomer	C ₂₅ H ₂₆ FN ₃ O	404.2	404.2
G7	ADB-5Br-INACA	C ₁₄ H ₁₇ BrN ₄ O ₂	353.1	353
G8	ADB-5'Br-BUTINACA	C ₁₈ H ₂₅ BrN ₄ O ₂	409.1	409
G9	CH-FUBIATA	C ₂₃ H ₂₅ FN ₂ O	365.2	365
G10	5-fluoro MPP-PICA phenylpropionic acid metabolite	C ₂₃ H ₂₅ FN ₂ O ₃	397.2	397.1
G11	CUMYL-INACA	C ₁₇ H ₁₇ N ₃ O	280.1	280.1
G12	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
H1	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>
H2	ADB-5'Br-DECINACA	C ₂₄ H ₃₇ BrN ₄ O ₂	493.2	493.2
H3	ADB-5'Br-PINACA	C ₁₉ H ₂₇ BrN ₄ O ₂	423.1	423.1
H4	ADB-5'Br-4en-PINACA	C ₁₉ H ₂₅ BrN ₄ O ₂	421.1	421.1
H5	ADB-INACA	C ₁₄ H ₁₈ N ₄ O ₂	275.2	275.2
H6	ADB-IATA	C ₁₆ H ₂₁ N ₃ O ₂	288.2	288.2
H7	CH-HEXIATA	C ₂₂ H ₃₂ N ₂ O	341.3	341.1
H8	CHM-FUBIATA	C ₂₄ H ₂₇ FN ₂ O	379.2	379.2
H9	5-fluoro MDMA-7-PAICA butanoic acid metabolite	C ₁₉ H ₂₆ FN ₃ O ₃	364.2	364
H10	4-cyano AB-BUTICA	C ₁₉ H ₂₄ N ₄ O ₂	341.2	341.2
H11	MN-18 N-(5-hydroxypentyl) metabolite	C ₂₃ H ₂₃ N ₃ O ₂	374.2	374.1
H12	<i>Unused</i>	<i>Unused</i>	<i>Unused</i>	<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

techserv@caymanchem.com
www.caymanchem.com