

CERTIFICATE of ANALYSIS



Tryptamine Analytical Standards Panel

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

Item No.:	33685
Batch No.:	0682557
Size:	100 µg/well
Expiry Date:	25SEP2026 (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 96-well Matrix plates containing individual tryptamine and tryptamine-class standards. The individual well locations for each unique standard are listed in the table below.

Traceability Statement

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

Quality Testing

A single kit was pulled randomly from the batch and high resolution accurate mass analysis was performed to verify the contents of each vial position. A summary of results is reported below as the average ppm mass error $\pm$ standard deviation determined for all analytes.

Test	Result	Specification	Meets Specification
Average mass error	3.08 $\pm$ 1.02 ppm	$\pm$ 10 ppm	Yes

Individual results are reported in the table on pages 2-5.

Approval: 

Title: Director of Analytical Chemistry

Certification Date: 25SEP2023

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)	Calculated [M+H] ⁺ m/z	Measured [M+H] ⁺ m/z	Mass Accuracy (ppm)
A1	Unused	Unused	Unused	Unused	Unused
A2	Tryptamine	C ₁₀ H ₁₂ N ₂	161.1079	161.1071	5.14
A3	α-methyl Tryptamine (hydrochloride)	C ₁₁ H ₁₄ N ₂	175.1235	175.1227	4.61
A4	N,N-DET (hydrochloride)	C ₁₄ H ₂₀ N ₂	217.1705	217.1696	4.27
A5	N-methyl Tryptamine	C ₁₁ H ₁₄ N ₂	175.1235	175.1227	4.90
A6	4-hydroxy MET	C ₁₃ H ₁₈ N ₂ O	219.1497	219.1488	4.21
A7	Gramine	C ₁₁ H ₁₄ N ₂	175.1235	175.1227	4.84
A8	5-methoxy DPT	C ₁₇ H ₂₆ N ₂ O	275.2123	275.2113	3.83
A9	7-fluoro Tryptamine (hydrochloride)	C ₁₀ H ₁₁ FN ₂	179.0985	179.0976	4.72
A10	DiPT	C ₁₆ H ₂₄ N ₂	245.2018	245.2007	4.36
A11	α-Ethyltryptamine	C ₁₂ H ₁₆ N ₂	189.1392	189.1383	4.85
A12	Unused	Unused	Unused	Unused	Unused
B1	Unused	Unused	Unused	Unused	Unused
B2	N,N-DMT (fumarate)	C ₁₂ H ₁₆ N ₂	189.1392	189.1382	5.01
B3	5-hydroxy-N-methyl Tryptamine (oxalate)	C ₁₁ H ₁₄ N ₂ O	191.1184	191.1174	5.46
B4	5-fluoro AMT (hydrochloride)	C ₁₁ H ₁₃ FN ₂	193.1141	193.1131	5.36
B5	4-hydroxy DET	C ₁₄ H ₂₀ N ₂ O	233.1654	233.1642	4.99
B6	5-chloro Tryptamine (hydrochloride)	C ₁₀ H ₁₁ ClN ₂	195.0689	195.0680	4.85
B7	4-methyl-α-Ethyltryptamine	C ₁₃ H ₁₈ N ₂	203.1548	203.1537	5.65
B8	5-methoxy MET	C ₁₄ H ₂₀ N ₂ O	233.1654	233.1643	4.56
B9	2-methyl DMT	C ₁₃ H ₁₈ N ₂	203.1548	203.1539	4.47
B10	5-methoxy EiPT	C ₁₆ H ₂₄ N ₂ O	261.1967	261.1955	4.57
B11	5-hydroxy DMT (hydrochloride)	C ₁₂ H ₁₆ N ₂ O	205.1341	205.1332	4.55
B12	Unused	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)	Calculated [M+H] ⁺ m/z	Measured [M+H] ⁺ m/z	Mass Accuracy (ppm)
C1	Unused	Unused	Unused	Unused	Unused
C2	4-hydroxy DMT	C ₁₂ H ₁₆ N ₂ O	205.1341	205.1330	5.13
C3	MiPT	C ₁₄ H ₂₀ N ₂	217.1705	217.1695	4.64
C4	5-IT	C ₁₁ H ₁₄ N ₂	175.1235	175.1227	4.90
C5	4-acetoxy MiPT (hydrochloride)	C ₁₆ H ₂₂ N ₂ O ₂	275.1760	275.1747	4.68
C6	6-IT	C ₁₁ H ₁₄ N ₂	175.1235	175.1226	5.13
C7	5-methoxy DMT	C ₁₃ H ₁₈ N ₂ O	219.1497	219.1487	4.80
C8	4-hydroxy DiPT (hydrochloride)	C ₁₆ H ₂₄ N ₂ O	261.1967	261.1955	4.49
C9	4-methoxy DMT	C ₁₃ H ₁₈ N ₂ O	219.1497	219.1487	4.67
C10	6-methoxy DiPT (hydrochloride)	C ₁₇ H ₂₆ N ₂ O	275.2123	275.2112	4.33
C11	5,7-dichloro Tryptamine (hydrochloride)	C ₁₀ H ₁₀ Cl ₂ N ₂	229.0299	229.0288	4.91
C12	Unused	Unused	Unused	Unused	Unused
D1	Unused	Unused	Unused	Unused	Unused
D2	4-hydroxy McPT (hydrochloride)	C ₁₄ H ₁₈ N ₂ O	231.1497	231.1488	4.12
D3	Norpsilocin	C ₁₁ H ₁₄ N ₂ O	191.1184	191.1176	4.57
D4	EPT	C ₁₅ H ₂₂ N ₂	231.1861	231.1851	4.49
D5	4-hydroxy MiPT	C ₁₄ H ₂₀ N ₂ O	233.1654	233.1643	4.86
D6	Ibogaine (hydrochloride)	C ₂₀ H ₂₆ N ₂ O	311.2123	311.2112	3.61
D7	4-hydroxy MPT	C ₁₄ H ₂₀ N ₂ O	233.1654	233.1644	4.26
D8	MET (oxalate)	C ₁₃ H ₁₈ N ₂	203.1548	203.1539	4.47
D9	6-fluoro DET	C ₁₄ H ₁₉ FN ₂	235.1611	235.1601	4.07
D10	N,N-DMT	C ₁₂ H ₁₆ N ₂	189.1392	189.1384	4.32
D11	DPT (hydrochloride)	C ₁₆ H ₂₄ N ₂	245.2018	245.2008	4.11
D12	Unused	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)	Calculated [M+H] ⁺ m/z	Measured [M+H] ⁺ m/z	Mass Accuracy (ppm)
E1	Unused	Unused	Unused	Unused	Unused
E2	4-acetoxy DMT (hydrochloride)	C ₁₄ H ₁₈ N ₂ O ₂	247.1447	247.1438	3.59
E3	5-methoxy MiPT	C ₁₅ H ₂₂ N ₂ O	247.1810	247.1800	4.14
E4	MALT	C ₁₄ H ₁₈ N ₂	215.1548	215.1540	3.99
E5	4-methoxy MiPT (hydrochloride)	C ₁₅ H ₂₂ N ₂ O	247.1810	247.1801	3.94
E6	5-methoxy DiBT (hydrochloride)	C ₁₉ H ₃₀ N ₂ O	303.2436	303.2427	3.11
E7	4-hydroxy EPT (hemifumarate)	C ₁₅ H ₂₂ N ₂ O	247.1810	247.1801	3.69
E8	5-methoxy-α-Ethyltryptamine	C ₁₃ H ₁₈ N ₂ O	219.1497	219.1488	4.49
E9	5-methoxy EPT (hydrochloride)	C ₁₆ H ₂₄ N ₂ O	261.1967	261.1958	3.27
E10	5-methoxy AMT	C ₁₂ H ₁₆ N ₂ O	205.1341	205.1332	4.16
E11	4-hydroxy DPT	C ₁₆ H ₂₄ N ₂ O	261.1967	261.1957	3.84
E12	Unused	Unused	Unused	Unused	Unused
F1	Unused	Unused	Unused	Unused	Unused
F2	5-methoxy DALT	C ₁₇ H ₂₂ N ₂ O	271.1810	271.1799	4.21
F3	4-hydroxy MALT	C ₁₄ H ₁₈ N ₂ O	231.1497	231.1489	3.86
F4	4-acetoxy DET (hydrochloride)	C ₁₆ H ₂₂ N ₂ O ₂	275.1760	275.1748	4.35
F5	N-acetyl-5-Hydroxytryptamine	C ₁₂ H ₁₄ N ₂ O ₂	219.1134	219.1123	4.69
F6	PiPT	C ₁₆ H ₂₄ N ₂	245.2018	245.2009	3.74
F7	5-methoxy DiPT	C ₁₇ H ₂₆ N ₂ O	275.2123	275.2113	3.68
F8	5-Hydroxytryptamine (hydrochloride)	C ₁₀ H ₁₂ N ₂ O	177.1028	177.1020	4.42
F9	4-methoxy DiPT (hydrochloride)	C ₁₇ H ₂₆ N ₂ O	275.2123	275.2114	3.54
F10	5-chloro DMT (hydrochloride)	C ₁₂ H ₁₅ ClN ₂	223.1002	223.0992	4.60
F11	7-methoxy DiPT (hydrochloride)	C ₁₇ H ₂₆ N ₂ O	275.2123	275.2114	3.35
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 1 Continued

Well	Item Name	Molecular Formula (free analyte)	Calculated [M+H] ⁺ m/z	Measured [M+H] ⁺ m/z	Mass Accuracy (ppm)
G1	Unused	Unused	Unused	Unused	Unused
G2	Psilocybin	C ₁₂ H ₁₇ N ₂ O ₄ P	285.1004	285.0995	3.32
G3	4-acetoxy MET (hydrochloride)	C ₁₅ H ₂₀ N ₂ O ₂	261.1603	261.1593	3.74
G4	5-methoxy DET	C ₁₅ H ₂₂ N ₂ O	247.1810	247.1800	4.10
G5	5-methoxy DBT	C ₁₉ H ₃₀ N ₂ O	303.2436	303.2428	2.75
G6	Unused	Unused	Unused	Unused	Unused
G7	Unused	Unused	Unused	Unused	Unused
G8	Unused	Unused	Unused	Unused	Unused
G9	Unused	Unused	Unused	Unused	Unused
G10	Unused	Unused	Unused	Unused	Unused
G11	Unused	Unused	Unused	Unused	Unused
G12	Unused	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



Revision History

Revision No.	Date	Reason for Revision
01	25SEP2023	Initial version

Disclaimers

Material Safety Data

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some but not all of the information required for the safe and proper use of this material. Before use, review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

Warranty and Limitation of Remedy

Cayman Chemical Company makes no warranty or guarantee of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specification.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver of Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitations of Remedy located on our website and in our catalog.

This Certificate shall not be reproduced except in full, without written approval from the Cayman Chemical Company.

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