

PRODUCT INFORMATION



Phenethylamine Analytical Standards Panel

Item No. 31617

Storage: -20°C
Stability: ≥5 years

Laboratory Procedures

Uncap each vial to be used. Add 500 µl of methanol to each vial. This will provide a 200 µg/ml standard solution for each analyte. Re-cap the vials and place plate on a plate mixer or vortexer. Mix for a minimum of 15 minutes to ensure full reconstitution.

The plate included in this panel contains a residual amount of glycerol to aid in the reconstitution of the analyte in methanol. Recovery in methanol has been validated for all analytes on the plate. Recovery from other solvents has not been verified.

Description

The Phenethylamine Analytical Standards Panel contains 71 analytical reference materials and standards categorized as phenethylamines. These compounds are supplied at 100 µg/well in a 96-well Matrix™ tube rack format for rapid screening or building in-house spectral libraries. The plate included in this panel contains a residual amount of glycerol to aid in the reconstitution of the analyte in methanol. Recovery in methanol has been validated for all analytes on the plate. Recovery from other solvents has not been verified. Please review the product insert for a full list of targets. The Phenethylamine Analytical Standards Panel 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. Please review the product insert for a full list of targets.

Panel Contents

Plate	Well	Contents	Item Number	Molecular Formula	CAS Number
1	A1	Unused			
1	A2	2-Phenethylamine (hydrochloride)	14206	C ₈ H ₁₁ N • HCl	156-28-5
1	A3	25P-NBOMe (hydrochloride)	9001991	C ₂₁ H ₂₉ NO ₃ • HCl	1539266-43-7
1	A4	N,N-Dimethylphenethylamine (hydrochloride)	29754	C ₁₀ H ₁₅ N • HCl	10275-21-5
1	A5	5-AEDB (hydrochloride)	18351	C ₁₀ H ₁₃ NO • HCl	796869-33-5
1	A6	Hordenine	21626	C ₁₀ H ₁₅ NO	539-15-1
1	A7	TDIQ (hydrochloride)	16083	C ₁₀ H ₁₁ NO ₂ • HCl	15052-05-8
1	A8	Homarylamine (hydrochloride)	21351	C ₁₀ H ₁₃ NO ₂ • HCl	533-10-8
1	A9	25T7-NBOMe (hydrochloride)	9001988	C ₂₁ H ₂₉ NO ₃ S • HCl	1539266-55-1
1	A10	3,4-Dimethoxyphenethylamine (hydrochloride)	14141	C ₁₀ H ₁₅ NO ₂ • HCl	635-85-8
1	A11	N-methyl Homarylamine (hydrochloride)	22239	C ₁₁ H ₁₅ NO ₂ • HCl	136366-39-7
1	A12	Unused			
1	B1	Unused			
1	B2	2C-D (hydrochloride)	11888	C ₁₁ H ₁₇ NO ₂ • HCl	25505-65-1
1	B3	25B-NB4OMe (hydrochloride)	19499	C ₁₈ H ₂₂ BrNO ₃ • HCl	1566571-56-9
1	B4	2C-G (hydrochloride)	14064	C ₁₂ H ₁₉ NO ₂ • HCl	327175-14-4
1	B5	Mescaline (hydrochloride)	11950	C ₁₁ H ₁₇ NO ₃ • HCl	832-92-8
1	B6	2C-C (hydrochloride)	11735	C ₁₀ H ₁₄ ClNO ₂ • HCl	88441-15-0
1	B7	25G-NBOMe (hydrochloride)	14065	C ₂₀ H ₂₇ NO ₃ • HCl	1797132-54-7
1	B8	2C-iP (hydrochloride)	18281	C ₁₃ H ₂₁ NO ₂ • HCl	1498978-47-4
1	B9	Escaline (hydrochloride)	14107	C ₁₂ H ₁₉ NO ₃ • HCl	3166-82-3

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, 09/30/2024

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	Molecular Formula	CAS Number
1	B10	25H-NBOMe (hydrochloride)	12061	$C_{18}H_{23}NO_3 \cdot HCl$	1566571-52-5
1	B11	N-methyl Mescaline (hydrochloride)	30541	$C_{12}H_{19}NO_3 \cdot HCl$	6308-81-2
1	B12	<i>Unused</i>			
1	C1	<i>Unused</i>			
1	C2	(±)-Lefetamine (hydrochloride)	15566	$C_{16}H_{19}N \cdot HCl$	24301-90-4
1	C3	2C-N	11890	$C_{10}H_{14}N_2O_4$	261789-00-8
1	C4	2C-T (hydrochloride)	13957	$C_{11}H_{17}NO_2S \cdot HCl$	61638-10-6
1	C5	Allylescaline (hydrochloride)	9001982	$C_{13}H_{19}NO_3 \cdot HCl$	39201-76-8
1	C6	(S)-Desoxy-D2PM (hydrochloride)	9001095	$C_{17}H_{19}N \cdot HCl$	188398-87-0
1	C7	Proscaline (hydrochloride)	19926	$C_{13}H_{21}NO_3 \cdot HCl$	61367-69-9
1	C8	2C-T-2 (hydrochloride)	11891	$C_{12}H_{19}NO_2S \cdot HCl$	681160-71-4
1	C9	2C-TFM (hydrochloride)	14471	$C_{11}H_{14}F_3NO_2 \cdot HCl$	159277-13-1
1	C10	25I-NBOMe 3-methoxy isomer (hydrochloride)	14784	$C_{18}H_{22}INO_3 \cdot HCl$	1566571-63-8
1	C11	2C-T-4 (hydrochloride)	11892	$C_{13}H_{21}NO_2S \cdot HCl$	868738-44-7
1	C12	<i>Unused</i>			
1	D1	<i>Unused</i>			
1	D2	2C-B (hydrochloride)	11734	$C_{10}H_{14}BrNO_2 \cdot HCl$	56281-37-9
1	D3	2-bromo-4,5-Dimethoxyphenethylamine (hydrochloride)	29386	$C_{10}H_{14}BrNO_2 \cdot HCl$	103275-23-6
1	D4	Sibutramine (hydrochloride)	14476	$C_{17}H_{26}NCl \cdot HCl$	84485-00-7
1	D5	2C-B-FLY (hydrochloride)	11473	$C_{12}H_{14}BrNO_2 \cdot HCl$	178557-21-6
1	D6	PRL-8-53 (hydrochloride)	26208	$C_{18}H_{21}NO_2 \cdot HCl$	51352-87-5
1	D7	25H-NBOH (hydrochloride)	22033	$C_{17}H_{21}NO_3 \cdot HCl$	2595087-21-9
1	D8	25H-NBOMe imine analog	12062	$C_{18}H_{21}NO_3$	1566571-78-5
1	D9	BD 1008 (hydrobromide)	9001336	$C_{15}H_{22}Cl_2N_2 \cdot 2HBr$	138356-09-9
1	D10	Metaescaline (hydrochloride)	19927	$C_{12}H_{19}NO_3 \cdot HCl$	90132-32-4
1	D11	25H-NB4OMe (hydrochloride)	19365	$C_{18}H_{23}NO_3 \cdot HCl$	1566571-54-7
1	D12	<i>Unused</i>			
1	E1	<i>Unused</i>			
1	E2	2C-I (hydrochloride)	11113	$C_{10}H_{14}INO_2 \cdot HCl$	64584-32-3
1	E3	25D-NBOMe (hydrochloride)	11946	$C_{19}H_{25}NO_3 \cdot HCl$	1539266-35-7
1	E4	25C-NBOH (hydrochloride)	14815	$C_{17}H_{20}ClNO_3 \cdot HCl$	1539266-20-0
1	E5	25C-NBF (hydrochloride)	15969	$C_{17}H_{19}ClFNO_2 \cdot HCl$	1539266-21-1
1	E6	RH-34	14445	$C_{18}H_{19}N_3O_3$	1028307-48-3
1	E7	2C-P (hydrochloride)	11696	$C_{13}H_{21}NO_2 \cdot HCl$	1359704-27-0
1	E8	25E-NBOMe (hydrochloride)	14727	$C_{20}H_{27}NO_3 \cdot HCl$	1539266-39-1
1	E9	Mescaline NBOMe (hydrochloride)	9001990	$C_{19}H_{25}NO_4 \cdot HCl$	2748304-49-4
1	E10	25C-NB3OMe (hydrochloride)	18560	$C_{18}H_{22}ClNO_3 \cdot HCl$	1566571-57-0
1	E11	25I-NBOMe (hydrochloride)	9001128	$C_{18}H_{22}INO_3 \cdot HCl$	1043868-97-8
1	E12	<i>Unused</i>			
1	F1	<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	Molecular Formula	CAS Number
1	F2	25iP-NBOMe (hydrochloride)	18170	C ₂₁ H ₂₉ NO ₃ • HCl	2725318-60-3
1	F3	N-methyl-2-Phenylpropan-1-amine (hydrochloride)	18563	C ₁₀ H ₁₅ N • HCl	5969-39-1
1	F4	25N-NBOMe (hydrochloride)	14911	C ₁₈ H ₂₂ N ₂ O ₅ • HCl	1566571-65-0
1	F5	25T-NBOMe (hydrochloride)	14562	C ₁₉ H ₂₅ NO ₃ S • HCl	1539266-47-1
1	F6	25T2-NBOMe (hydrochloride)	9002690	C ₂₀ H ₂₇ NO ₃ S • HCl	1539266-51-7
1	F7	25B-NBOH (hydrochloride)	18328	C ₁₇ H ₂₀ BrNO ₃ • HCl	1539266-16-4
1	F8	25B-NBF (hydrochloride)	15967	C ₁₇ H ₁₉ BrFNO ₂ • HCl	1539266-17-5
1	F9	2C-H (hydrochloride)	11889	C ₁₀ H ₁₅ NO ₂ • HCl	3166-74-3
1	F10	25T4-NBOMe (hydrochloride)	9001989	C ₂₁ H ₂₉ NO ₃ S • HCl	1566571-73-0
1	F11	25B-NBOMe (hydrochloride)	12047	C ₁₈ H ₂₂ BrNO ₃ • HCl	1539266-15-3
1	F12	<i>Unused</i>			
1	G1	<i>Unused</i>			
1	G2	25B-NB3OMe (hydrochloride)	19364	C ₁₈ H ₂₂ BrNO ₃ • HCl	1566571-55-8
1	G3	2C-E (hydrochloride)	10395	C ₁₂ H ₁₉ NO ₂ • HCl	923013-67-6
1	G4	30C-NBOMe (hydrochloride)	14528	C ₂₀ H ₂₆ ClNO ₅ • HCl	2749391-65-7
1	G5	25I-NBOH (hydrochloride)	14909	C ₁₇ H ₂₀ INO ₃ • HCl	1539266-12-0
1	G6	25I-NBF (hydrochloride)	14066	C ₁₇ H ₁₉ FINO ₂ • HCl	1539266-13-1
1	G7	25I-NBOMe imine analog	11968	C ₁₈ H ₂₀ INO ₃	2177287-33-9
1	G8	2C-T-7 (hydrochloride)	11200	C ₁₃ H ₂₁ NO ₂ S • HCl	850140-15-7
1	G9	25I-NBOMe 4-methoxy isomer (hydrochloride)	14785	C ₁₈ H ₂₂ INO ₃ • HCl	1566571-64-9
1	G10	25C-NBOMe (hydrochloride)	9001096	C ₁₈ H ₂₂ ClNO ₃ • HCl	1539266-19-7
1	G11	25I-NBMD (hydrochloride)	11947	C ₁₈ H ₂₀ INO ₄ • HCl	1539266-14-2
1	G12	<i>Unused</i>			
1	H1	<i>Unused</i>			
1	H2	N-acetyl 25I-NBOMe	24806	C ₂₀ H ₂₄ INO ₄	2748301-01-9

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