

PRODUCT INFORMATION



Atypical Stimulants Analytical Standards Panel

Item No. 31618

Storage: -20°C
Stability: ≥3 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 Atypical Stimulants Analytical Standards Panel contains 126 analytical reference materials and standards categorized as atypical stimulants, including arylcyclohexylamines, piperazines, aminoindanes, opioids, nootropics, and precursors in the synthesis of stimulants. 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 Atypical Stimulants Analytical Standards Panel contains compounds regulated as Schedule I-IV 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	Molecular Formula	CAS Number
1	A1	<i>Unused</i>			
1	A2	Piperazine	24019	C ₄ H ₁₀ N ₂	110-85-0
1	A3	1,3-Dimethylbutylamine (hydrochloride)	21928	C ₆ H ₁₅ N • HCl	71776-70-0
1	A4	Octodrine (hydrochloride)	21927	C ₈ H ₁₉ N • HCl	5984-59-8
1	A5	2-AI (hydrochloride)	11048	C ₉ H ₁₁ N • HCl	2338-18-3
1	A6	Phenylacetone	16103	C ₉ H ₁₀ O	103-79-7
1	A7	Phenylacetic Acid	18709	C ₈ H ₈ O ₂	103-82-2
1	A8	N-methyl-2-AI (hydrochloride)	14897	C ₁₀ H ₁₃ N • HCl	10408-85-2
1	A9	Cathine	13868	C ₉ H ₁₃ NO	492-39-7
1	A10	1-(4-Methoxyphenyl)piperazine (hydrochloride)	14020	C ₁₁ H ₁₆ N ₂ O • 2HCl	38869-47-5
1	A11	α-Phenylacetoacetonitrile	21105	C ₁₀ H ₉ NO	4468-48-8
1	A12	<i>Unused</i>			
1	B1	<i>Unused</i>			
1	B2	Phenylpiperazine (hydrochloride)	11203	C ₁₀ H ₁₄ N ₂ • 2HCl	4004-95-9
1	B3	Phenyl-2-nitropropene	22185	C ₉ H ₉ NO ₂	705-60-2
1	B4	MEAI (hydrochloride)	18286	C ₁₀ H ₁₃ NO • HCl	81593-54-6
1	B5	1-Benzyl-2-methylpiperazine (hydrochloride)	23708	C ₁₂ H ₁₈ N ₂ • 2HCl	93138-53-5
1	B6	(1S,2S)-(+)-Pseudoephedrine	14209	C ₁₀ H ₁₅ NO	90-82-4
1	B7	3',4'-Dihydroxyphenylacetone	19522	C ₉ H ₁₀ O ₃	2503-44-8
1	B8	5-APDI (hydrochloride)	13899	C ₁₂ H ₁₇ N • HCl	152623-95-5
1	B9	BZP (hydrochloride)	11202	C ₁₁ H ₁₆ N ₂ • 2HCl	5321-63-1

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, 08/26/2024

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	BMK Glycidic Acid (sodium salt)	19683	C ₁₀ H ₉ O ₃ • Na	5449-12-7
1	B11	MDAI (hydrochloride)	9001102	C ₁₀ H ₁₁ NO ₂ • HCl	155344-90-4
1	B12	<i>Unused</i>			
1	C1	<i>Unused</i>			
1	C2	4,5-MDAI (hydrochloride)	9002653	C ₁₀ H ₁₁ NO ₂ • HCl	124399-90-2
1	C3	MMAI (hydrochloride)	13895	C ₁₁ H ₁₅ NO • HCl	132980-17-7
1	C4	Piperonyl methyl ketone	13892	C ₁₀ H ₁₀ O ₃	4676-39-5
1	C5	1-(4-Trifluoromethylphenyl)piperazine (hydrochloride)	21949	C ₁₁ H ₁₃ F ₃ N ₂ • 2HCl	2640680-83-5
1	C6	1-(4-Fluorophenyl)piperazine (hydrochloride)	11204	C ₁₀ H ₁₃ FN ₂ • 2HCl	64090-19-3
1	C7	1-(4-Chlorophenyl)piperazine	14481	C ₁₀ H ₁₃ ClN ₂	38212-33-8
1	C8	1-(2-Fluorophenyl)piperazine (hydrochloride)	21954	C ₁₀ H ₁₃ FN ₂ • 2HCl	76835-09-1
1	C9	(+)-Chloropseudoephedrine (hydrochloride)	9002347	C ₁₀ H ₁₄ ClN • HCl	94133-42-3
1	C10	Ecgonine (hydrochloride)	23314	C ₉ H ₁₅ NO ₃ • HCl	5796-31-6
1	C11	Deschloronorketamine (hydrochloride)	26401	C ₁₂ H ₁₅ NO • HCl	7015-20-5
1	C12	<i>Unused</i>			
1	D1	<i>Unused</i>			
1	D2	4,4'-Dimethylaminorex	15281	C ₁₁ H ₁₄ N ₂ O	1445569-01-6
1	D3	4-fluoro Deschloroketamine (hydrochloride)	29755	C ₁₃ H ₁₆ FNO • HCl	
1	D4	MBZP (hydrochloride)	11736	C ₁₂ H ₁₈ N ₂ • HCl	374898-00-7
1	D5	(+)-Ephedrine (hydrochloride)	14143	C ₁₀ H ₁₅ NO • HCl	24221-86-1
1	D6	1-Benzyl-3-methylpiperazine	23709	C ₁₂ H ₁₈ N ₂	3138-90-7
1	D7	Cocaine	16186	C ₁₇ H ₂₁ NO ₄	50-36-2
1	D8	Methyl 2-phenylacetoacetate	28352	C ₁₁ H ₁₂ O ₃	16648-44-5
1	D9	4-methoxy PCE (hydrochloride)	9001356	C ₁₅ H ₂₃ NO • HCl	1933-15-9
1	D10	DL-Norephedrine (hydrochloride)	14212	C ₉ H ₁₃ NO • HCl	154-41-6
1	D11	1-(3-Methoxyphenyl)piperazine	19273	C ₁₁ H ₁₆ N ₂ O	16015-71-7
1	D12	<i>Unused</i>			
1	E1	<i>Unused</i>			
1	E2	1-Piperidinocyclohexanecarbonitrile	20367	C ₁₂ H ₂₀ N ₂	3867-15-0
1	E3	Caffeine	22975	C ₈ H ₁₀ N ₄ O ₂	58-08-2
1	E4	4-fluoro BZP (hydrochloride)	11112	C ₁₁ H ₁₅ FN ₂ • 2HCl	199672-06-5
1	E5	3-Fluorophenmetrazine (hydrochloride)	17208	C ₁₁ H ₁₄ FNO • HCl	1803562-83-5
1	E6	1-(3-Chlorophenyl)piperazine (hydrochloride)	13949	C ₁₀ H ₁₃ ClN ₂ • HCl	13078-15-4
1	E7	1-(3-Fluorophenyl)piperazine (hydrochloride)	21953	C ₁₀ H ₁₃ FN ₂ • 2HCl	76835-10-4
1	E8	Deschloroketamine (hydrochloride)	18683	C ₁₃ H ₁₇ NO • HCl	4631-27-0
1	E9	BMK ethyl glycidate	9002713	C ₁₂ H ₁₄ O ₃	41232-97-7
1	E10	2-fluoro Deschloronorketamine (hydrochloride)	29118	C ₁₂ H ₁₄ FNO • HCl	1956327-93-7
1	E11	1-(2-Chloro-3-fluorophenyl)piperazine (hydrochloride)	29116	C ₁₀ H ₁₂ ClFN ₂ • HCl	
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	1-(3-Chloro-4-fluorophenyl)piperazine (hydrochloride)	30230	C ₁₀ H ₁₂ ClFN ₂ • HCl	95884-48-3
1	F3	Harmaline	10995	C ₁₃ H ₁₄ N ₂ O	304-21-2
1	F4	deschloro-N-ethyl-Ketamine (hydrochloride)	14105	C ₁₄ H ₁₉ NO • HCl	4551-92-2
1	F5	PCPr (hydrochloride)	9001362	C ₁₅ H ₂₃ N • HCl	1934-55-0
1	F6	Normethoxetamine (hydrochloride)	30757	C ₁₃ H ₁₇ NO ₂ • HCl	
1	F7	1-(3-Methylbenzyl)piperazine	24018	C ₁₂ H ₁₈ N ₂	5321-48-2
1	F8	Methyl 2-phenyl-2-(pyrrolidin-1-yl)acetate (hydrochloride)	30228	C ₁₃ H ₁₇ NO ₂ • HCl	27594-60-1
1	F9	3-hydroxy PCE (hydrochloride)	26877	C ₁₄ H ₂₁ NO • HCl	
1	F10	MDBP (hydrochloride)	19215	C ₁₂ H ₁₆ N ₂ O ₂ • 2HCl	38063-96-6
1	F11	PMK Glycidic Acid (sodium salt)	19685	C ₁₁ H ₉ O ₅ • Na	
1	F12	Unused			
1	G1	Unused			
1	G2	2-fluoro Deschloroketamine (hydrochloride)	19786	C ₁₃ H ₁₆ FNO • HCl	111982-49-1
1	G3	Benzoylpiperazine	24016	C ₁₁ H ₁₄ N ₂ O	13754-38-6
1	G4	3-fluoro Deschloroketamine (hydrochloride)	29756	C ₁₃ H ₁₆ FNO • HCl	2657761-24-3
1	G5	Norketamine (hydrochloride)	15787	C ₁₂ H ₁₄ ClNO • HCl	79499-59-5
1	G6	Tiletamine (hydrochloride)	19794	C ₁₂ H ₁₇ NOS • HCl	14176-50-2
1	G7	Theacrine	26855	C ₉ H ₁₂ N ₄ O ₃	2309-49-1
1	G8	Ephedrine (hydrochloride)	18327	C ₁₆ H ₁₉ N • HCl	6272-97-5
1	G9	Rolicyclidine (hydrochloride)	19263	C ₁₆ H ₂₃ N • HCl	1934-48-1
1	G10	2,3-Dichlorophenylpiperazine (hydrochloride)	14296	C ₁₀ H ₁₂ Cl ₂ N ₂ • HCl	119532-26-2
1	G11	1-(3-Trifluoromethylphenyl)piperazine (hydrochloride)	11205	C ₁₁ H ₁₃ F ₃ N ₂ • 2HCl	76835-14-8
1	G12	Unused			
1	H1	Unused			
1	H2	1-(2-Trifluoromethylphenyl)piperazine (hydrochloride)	21948	C ₁₁ H ₁₃ F ₃ N ₂ • HCl	40160-26-7
1	H3	BMK Glycidic Acid	19682	C ₁₀ H ₁₀ O ₃	25547-51-7
1	H4	2-(3-(Trifluoromethyl)phenyl)piperazine (hydrochloride)	24420	C ₁₁ H ₁₃ F ₃ N ₂ • 2HCl	2517636-54-1
1	H5	2-methoxy Ketamine (hydrochloride)	12048	C ₁₄ H ₁₉ NO ₂ • HCl	6728-62-7
1	H6	(R)-Ketamine (hydrochloride)	16519	C ₁₃ H ₁₆ ClNO • HCl	33795-24-3
1	H7	Methoxmetamine (hydrochloride)	17168	C ₁₄ H ₁₉ NO ₂ • HCl	1781829-56-8
1	H8	2'-hydroxy Cocaine	20062	C ₁₇ H ₂₁ NO ₅	89339-17-3
1	H9	(S)-(+)-MT-45 (hydrochloride)	15128	C ₂₄ H ₃₂ N ₂ • 2HCl	52694-54-9
1	H10	PMK methyl glycidate	21833	C ₁₂ H ₁₂ O ₅	13605-48-6
1	H11	1-(2,5-Dimethoxybenzyl)piperazine (hydrochloride)	24017	C ₁₃ H ₂₀ N ₂ O ₂ • HCl	374897-99-1
1	H12	Unused			
2	A1	Unused			
2	A2	Ketamine (hydrochloride)	11630	C ₁₃ H ₁₆ ClNO • HCl	1867-66-9
2	A3	(±)-threo-Methylphenidate (hydrochloride)	11639	C ₁₄ H ₁₉ NO ₂ • HCl	23655-65-4
2	A4	(S)-Ketamine (hydrochloride)	9001961	C ₁₃ H ₁₆ ClNO • HCl	33643-47-9
2	A5	(2R,6R)-Hydroxynorketamine (hydrochloride)	19603	C ₁₂ H ₁₄ ClNO ₂ • HCl	1430202-69-9

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

PRODUCT INFORMATION



Plate	Well	Contents	Item Number	Molecular Formula	CAS Number
2	A6	1-(4-Bromophenyl)piperazine (hydrochloride)	30229	C ₁₀ H ₁₃ BrN ₂ • HCl	68104-62-1
2	A7	PCP (hydrochloride)	14276	C ₁₇ H ₂₅ N • HCl	956-90-1
2	A8	Norcocaine (hydrochloride)	28492	C ₁₆ H ₁₉ NO ₄ • HCl	61585-22-6
2	A9	(±)- <i>threo</i> -Ethylphenidate (hydrochloride)	13896	C ₁₅ H ₂₁ NO ₂ • HCl	214149-46-9
2	A10	(±)- <i>threo</i> -Isopropylphenidate (hydrochloride)	19264	C ₁₆ H ₂₃ NO ₂ • HCl	1262795-94-7
2	A11	PCEEA (hydrochloride)	9001365	C ₁₆ H ₂₅ NO • HCl	1798021-89-2
2	A12	Unused			
2	B1	Unused			
2	B2	Tenocyclidine (hydrochloride)	17014	C ₁₅ H ₂₃ NS • HCl	1867-65-8
2	B3	PMK ethyl glycidate	21990	C ₁₃ H ₁₄ O ₅	28578-16-7
2	B4	N-Ethylnorketamine (hydrochloride)	18591	C ₁₄ H ₁₈ CINO • HCl	2525095-53-6
2	B5	(±)- <i>threo</i> -4-Fluoromethylphenidate (hydrochloride)	9002651	C ₁₄ H ₁₈ FNO ₂ • HCl	2703832-68-0
2	B6	2-DPMP (hydrochloride)	11481	C ₁₈ H ₂₁ N • HCl	5807-81-8
2	B7	D2PM (hydrochloride)	11160	C ₁₇ H ₁₉ NO • HCl	172152-19-1
2	B8	Dextrorphan (tartrate)	15886	C ₁₇ H ₂₃ NO • C ₄ H ₆ O ₆	143-98-6
2	B9	5-IAI (hydrochloride)	11035	C ₉ H ₁₀ IN • HCl	1782044-60-3
2	B10	3-methoxy PCP (hydrochloride)	11137	C ₁₈ H ₂₇ NO • HCl	91164-58-8
2	B11	1-(1-Phenylcyclohexyl)-4-hydroxypiperidine	20563	C ₁₇ H ₂₅ NO	60232-85-1
2	B12	Unused			
2	C1	Unused			
2	C2	<i>trans</i> -4-phenyl-4-Piperidinocyclohexanol	20565	C ₁₇ H ₂₅ NO	78165-07-8
2	C3	Methoxpropamine (hydrochloride)	30248	C ₁₆ H ₂₃ NO ₂ • HCl	
2	C4	PCMPA (hydrochloride)	9001363	C ₁₆ H ₂₅ NO • HCl	1934-63-0
2	C5	(±)- <i>threo</i> -Propylphenidate (hydrochloride)	19268	C ₁₆ H ₂₃ NO ₂ • HCl	2514950-54-8
2	C6	3- <i>p</i> -FBT (hydrochloride)	18057	C ₁₅ H ₁₈ FNO ₂ • HCl	
2	C7	N-Propylnorketamine (hydrochloride)	20272	C ₁₅ H ₂₀ CINO • HCl	2749326-65-4
2	C8	(±)- <i>threo</i> -4-Fluoroethylphenidate (hydrochloride)	20020	C ₁₅ H ₂₀ FNO ₂ • HCl	2740569-99-5
2	C9	Trimetazidine (hydrochloride)	18165	C ₁₄ H ₂₂ N ₂ O ₃ • 2HCl	13171-25-0
2	C10	1,4-Dibenzylpiperazine (hydrochloride)	11206	C ₁₈ H ₂₂ N ₂ • 2HCl	2298-55-7
2	C11	Pipradrol (hydrochloride)	14783	C ₁₈ H ₂₁ NO • HCl	71-78-3
2	C12	Unused			
2	D1	Unused			
2	D2	4-methoxy PCP (hydrochloride)	11136	C ₁₈ H ₂₇ NO • HCl	2185-93-5
2	D3	2C-B-BZP (hydrochloride)	14793	C ₁₃ H ₁₉ BrN ₂ O ₂ • 2HCl	2748622-68-4
2	D4	3-hydroxy PCP (hydrochloride)	14421	C ₁₇ H ₂₅ NO • HCl	79295-51-5
2	D5	2-bromo Deschloroketamine	30367	C ₁₃ H ₁₆ BrNO	120807-70-7
2	D6	(±)- <i>threo</i> -Methylnaphthidate (hydrochloride)	17970	C ₁₈ H ₂₁ NO ₂ • HCl	219915-69-2
2	D7	Diclofensine (hydrochloride)	15355	C ₁₇ H ₁₇ Cl ₂ NO • HCl	34041-84-4
2	D8	Adrafinil	19773	C ₁₅ H ₁₅ NO ₃ S	63547-13-7
2	D9	Methoxetamine (hydrochloride)	11139	C ₁₅ H ₂₁ NO ₂ • HCl	1239908-48-5
2	D10	Benzoylcegonine	9001495	C ₁₆ H ₁₉ NO ₄	519-09-5

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
2	D11	RTI-31	17977	C ₁₆ H ₂₀ ClNO ₂	130342-80-2
2	D12	Unused			
2	E1	Unused			
2	E2	(±)-threo-Ethyl-naphthidate (hydrochloride)	17954	C ₁₉ H ₂₃ NO ₂ • HCl	
2	E3	Benocyclidine	11737	C ₁₉ H ₂₅ NS	112726-66-6
2	E4	(±)-threo-3,4-Dichloromethylphenidate (hydrochloride)	18588	C ₁₄ H ₁₇ Cl ₂ NO ₂ • HCl	214149-42-5
2	E5	Scopolamine (hydrobromide)	14108	C ₁₇ H ₂₁ NO ₄ • HBr	114-49-8
2	E6	MT-45 (hydrochloride)	14082	C ₂₄ H ₃₂ N ₂ • 2HCl	57314-55-3
2	E7	Bromantane	24688	C ₁₆ H ₂₀ BrN	87913-26-6
2	E8	Unused			
2	E9	Unused			
2	E10	Unused			
2	E11	Unused			
2	E12	Unused			
2	F1	Unused			
2	F2	Unused			
2	F3	Unused			
2	F4	Unused			
2	F5	Unused			
2	F6	Unused			
2	F7	Unused			
2	F8	Unused			
2	F9	Unused			
2	F10	Unused			
2	F11	Unused			
2	F12	Unused			
2	G1	Unused			
2	G2	Unused			
2	G3	Unused			
2	G4	Unused			
2	G5	Unused			
2	G6	Unused			
2	G7	Unused			
2	G8	Unused			
2	G9	Unused			
2	G10	Unused			
2	G11	Unused			
2	G12	Unused			
2	H1	Unused			
2	H2	Unused			
2	H3	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	Molecular Formula	CAS Number
2	H4	<i>Unused</i>			
2	H5	<i>Unused</i>			
2	H6	<i>Unused</i>			
2	H7	<i>Unused</i>			
2	H8	<i>Unused</i>			
2	H9	<i>Unused</i>			
2	H10	<i>Unused</i>			
2	H11	<i>Unused</i>			
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