

PRODUCT INFORMATION

Benzodiazepine Analytical Standards Panel

Item No. 26260

Storage: -20°C
Stability: ≥4 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 Benzodiazepine Analytical Standards Panel contains 53 analytical reference materials and standards categorized as benzodiazepines and benzodiazepine metabolites. 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.

Please review the product insert for a full list of targets. The Benzodiazepine Analytical Standards Panel contains compounds regulated as Schedule IV compounds in the United States and is regulated as a Schedule IV 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	α-hydroxy Midazolam	10385	C ₁₈ H ₁₃ ClFN ₃ O	59468-90-5
1	A3	Ketazolam	11449	C ₂₀ H ₁₇ ClN ₂ O ₃	27223-35-4
1	A4	3-hydroxy Phenazepam	11472	C ₁₅ H ₁₀ BrClN ₂ O ₂	70030-11-4
1	A5	Etizolam	12063	C ₁₇ H ₁₅ ClN ₄ S	40054-69-1
1	A6	Alprazolam	14255	C ₁₇ H ₁₃ ClN ₄	28981-97-7
1	A7	7-Aminoflunitrazepam	14258	C ₁₆ H ₁₄ FN ₃ O	34084-50-9
1	A8	Bromazepam	14260	C ₁₄ H ₁₀ BrN ₃ O	1812-30-2
1	A9	Clonazepam	14263	C ₁₅ H ₁₀ ClN ₃ O ₃	1622-61-3
1	A10	Pyrazolam	14901	C ₁₆ H ₁₂ BrN ₅	39243-02-2
1	A11	Flubromazepam	15157	C ₁₅ H ₁₀ BrFN ₂ O	2647-50-9
1	A12	<i>Unused</i>			
1	B1	<i>Unused</i>			
1	B2	α-hydroxy Alprazolam	15287	C ₁₇ H ₁₃ ClN ₄ O	37115-43-8
1	B3	Diclazepam	15554	C ₁₆ H ₁₂ Cl ₂ N ₂ O	2894-68-0
1	B4	4-hydroxy Alprazolam	15878	C ₁₇ H ₁₃ ClN ₄ O	30896-57-2
1	B5	Diazepam	15887	C ₁₆ H ₁₃ ClN ₂ O	439-14-5
1	B6	Estazolam	15889	C ₁₆ H ₁₁ ClN ₄	29975-16-4
1	B7	Lorazepam	15891	C ₁₅ H ₁₀ Cl ₂ N ₂ O ₂	846-49-1
1	B8	Nordiazepam	15898	C ₁₅ H ₁₁ ClN ₂ O	1088-11-5

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, 07/25/2024

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



Page 2 of 2

1	B9	Oxazepam	15909	C ₁₅ H ₁₁ ClN ₂ O ₂	604-75-1
1	B10	Temazepam	15918	C ₁₆ H ₁₃ ClN ₂ O ₂	846-50-4
1	B11	Delorazepam	15938	C ₁₅ H ₁₀ C ₂ N ₂ O	2894-67-9
1	B12	<i>Unused</i>			
1	C1	<i>Unused</i>			
1	C2	Flunitrazepam	16188	C ₁₆ H ₁₂ FN ₃ O ₃	1622-62-4
1	C3	Flurazepam	16190	C ₂₁ H ₂₃ ClFN ₃ O	17617-23-1
1	C4	Midazolam	16193	C ₁₈ H ₁₃ ClFN ₃	59467-70-8
1	C5	Nitrazepam	16196	C ₁₅ H ₁₁ N ₃ O ₃	146-22-5
1	C6	Flubromazolam	16435	C ₁₇ H ₁₂ BrFN ₄	612526-40-6
1	C7	Clonazolam	18173	C ₁₇ H ₁₂ ClN ₅ O ₂	33887-02-4
1	C8	Adinazolam	18205	C ₁₉ H ₁₈ ClN ₅	37115-32-5
1	C9	Halazepam	18290	C ₁₇ H ₁₂ ClF ₃ N ₂ O	23092-17-3
1	C10	Nimetazepam	18352	C ₁₆ H ₁₃ N ₃ O ₃	2011-67-8
1	C11	7-Aminonimetazepam	18353	C ₁₆ H ₁₅ N ₃ O	4959-16-4
1	C12	<i>Unused</i>			
1	D1	<i>Unused</i>			
1	D2	Prazepam	18358	C ₁₉ H ₁₇ ClN ₂ O	2955-38-6
1	D3	Desalkylflurazepam	18484	C ₁₅ H ₁₀ ClFN ₂ O	2886-65-9
1	D4	N-Desmethyleclonazepam	18564	C ₁₅ H ₁₁ ClN ₂ O ₂	22316-55-8
1	D5	Cinolazepam	20529	C ₁₈ H ₁₃ ClFN ₃ O ₂	75696-02-5
1	D6	Fludiazepam	22424	C ₁₆ H ₁₂ ClFN ₂ O	3900-31-0
1	D7	Triazolam	22577	C ₁₇ H ₁₂ Cl ₂ N ₄	28911-01-5
1	D8	Nitrazolam	22620	C ₁₇ H ₁₃ N ₅ O ₂	28910-99-8
1	D9	Bromazolam	22665	C ₁₇ H ₁₃ BrN ₄	71368-80-4
1	D10	Methylclonazepam	22747	C ₁₆ H ₁₂ ClN ₃ O ₃	5527-71-9
1	D11	Cloniprazepam	23106	C ₁₉ H ₁₆ ClN ₃ O ₃	1998158-84-1
1	D12	<i>Unused</i>			
1	E1	<i>Unused</i>			
1	E2	Flunitrazolam	23108	C ₁₇ H ₁₂ FN ₅ O ₂	2243815-18-9
1	E3	Difludiazepam	23507	C ₁₆ H ₁₁ ClF ₂ N ₂ O	39080-67-6
1	E4	Fluclozepam	23524	C ₁₅ H ₁₀ ClFN ₄ S	54123-15-8
1	E5	1-demethyl Phenazolam	23525	C ₁₆ H ₁₀ BrClN ₄	919973-69-6
1	E6	Flutoprazepam	23543	C ₁₉ H ₁₆ ClFN ₂ O	25967-29-7
1	E7	Ethyl Loflazepate	23630	C ₁₈ H ₁₄ ClFN ₂ O ₃	29177-84-2
1	E8	Meclonazepam	25110	C ₁₆ H ₁₂ ClN ₃ O ₃	58662-84-3
1	E9	Flualprazolam	24481	C ₁₇ H ₁₂ ClFN ₄	28910-91-0
1	E10	Phenazolam	26700	C ₁₇ H ₁₂ BrClN ₄	87213-50-1
1	E11	8-Aminoclonazolam	27247	C ₁₇ H ₁₄ ClN ₅	71368-66-6
1	E12	<i>Unused</i>			
1	F1	<i>Unused</i>			
1	F2	Desmethyleclonazepam	27838	C ₁₅ H ₁₃ ClN ₂ OS	33671-37-3
1	F3	Phenazepam	9000849	C ₁₅ H ₁₀ BrClN ₂ O	51753-57-2
1	F4	N-Desmethyleflunitrazepam	9002548	C ₁₅ H ₁₀ FN ₃ O ₃	2558-30-7

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