

PRODUCT INFORMATION



Benzodiazepine Analytical Standards Panel

Item No. 26260

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 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, 04/10/2024

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-Desmethyloclobazam | 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-Amino-clonazolam | 27247 | C ₁₇ H ₁₄ ClN ₅ | 71368-66-6 |
| 1 | E12 | <i>Unused</i> | | | |
| 1 | F1 | <i>Unused</i> | | | |
| 1 | F2 | Desmethyloclozepam | 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-Desmethylflunitrazepam | 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