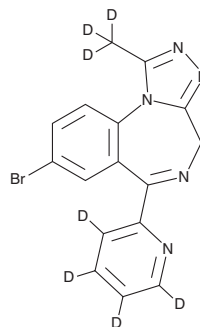


# PRODUCT INFORMATION



## Pyrazolam-d<sub>7</sub> Item No. 38176

**Formal Name:** 8-bromo-1-methyl-d<sub>3</sub>-6-(2-pyridinyl-d<sub>4</sub>)-4H-[1,2,4]triazolo[4,3-a][1,4]benzodiazepine  
**MF:** C<sub>16</sub>H<sub>5</sub>D<sub>7</sub>BrN<sub>5</sub>  
**FW:** 361.2  
**Chemical Purity:** ≥98% (Pyrazolam)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>7</sub>); ≤1% d<sub>0</sub>  
**UV/Vis.:** λ<sub>max</sub>: 223 nm  
**Supplied as:** A solution in acetonitrile  
**Storage:** -20°C  
**Stability:** ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description

Pyrazolam-d<sub>7</sub> (Item No. 38176) is intended for use as an internal standard for the quantification of pyrazolam (Item No. 14901) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Pyrazolam is categorized as a benzodiazepine.<sup>1</sup> Pyrazolam has been abused recreationally and has anxiolytic-like effects in mice.<sup>1,2</sup> This product is intended for research and forensic applications.

### References

1. Moosmann, B., Huppertz, L.M., Hutter, M., *et al.* Detection and identification of the designer benzodiazepine flubromazepam and preliminary data on its metabolism and pharmacokinetics. *J. Mass. Spectrom.* **48(11)**, 1150-1159 (2013).
2. Hester, J.B., Jr. and Von Voigtlander, P. 6-Aryl-4H-s-triazolo[4,3-a][1,4]benzodiazepines. Influence of 1-substitution on pharmacological activity. *J. Med. Chem.* **22(11)**, 1390-1398 (1979).

#### 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, 09/29/2023

#### 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