

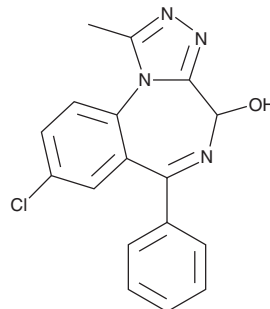
# PRODUCT INFORMATION



## 4-hydroxy Alprazolam

Item No. 15878

**CAS Registry No.:** 30896-57-2  
**Formal Name:** 8-chloro-1-methyl-6-phenyl-4H-[1,2,4]triazolo[4,3-a][1,4]benzodiazepin-4-ol  
**MF:** C<sub>17</sub>H<sub>13</sub>ClN<sub>4</sub>O  
**FW:** 324.8  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 223 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

Alprazolam is a benzodiazepine with anxiolytic and sedative actions.<sup>1</sup> 4-hydroxy Alprazolam is a major metabolite of alprazolam, produced by the actions of cytochrome P450 (CYP) isoforms CYP3A4 and CYP3A5.<sup>1-3</sup> It may be detected in serum or, in a glucuronidated form, in urine.<sup>2</sup> Alternatively, 4-hydroxyalprazolam can be further metabolized to benzophenone.<sup>2</sup>

### References

1. Park, J.-Y., Kim, K.-A., Park, P.-W., *et al.* Effect of CYP3A5\*3 genotype on the pharmacokinetics and pharmacodynamics of alprazolam in healthy subjects. *Clin. Pharmacol. Ther.* **79(6)**, 590-599 (2006).
2. Hirota, N., Ito, K., Iwatsubo, T., *et al.* *In vitro/in vivo* scaling of alprazolam metabolism by CYP3A4 and CYP3A5 in humans. *Biopharm. Drug Dispos.* **22(2)**, 53-71 (2001).
3. Williams, J.A., Ring, B.J., Cantrell, V.E., *et al.* Comparative metabolic capabilities of CYP3A4, CYP3A5, and CYP3A7. *Drug Metab. Dispos.* **30(8)**, 883-891 (2002).

#### 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, 06/28/2023

#### 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