

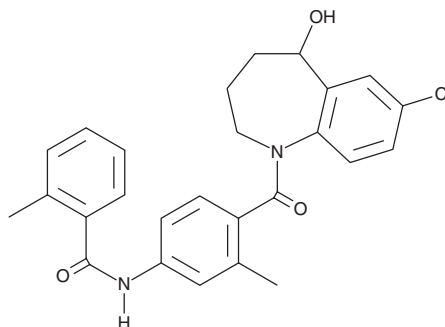
# PRODUCT INFORMATION



## Tolvaptan

Item No. 26301

**CAS Registry No.:** 150683-30-0  
**Formal Name:** N-[4-[(7-chloro-2,3,4,5-tetrahydro-5-hydroxy-1H-1-benzazepin-1-yl)carbonyl]-3-methylphenyl]-2-methyl-benzamide  
**Synonym:** OPC 41061  
**MF:** C<sub>26</sub>H<sub>25</sub>ClN<sub>2</sub>O<sub>3</sub>  
**FW:** 448.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 270 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

Tolvaptan (Item No. 26301) is an analytical reference standard categorized as a diuretic.<sup>1</sup> Formulations containing diuretics have been abused as performance-enhancing drugs and masking agents in sports doping.<sup>1,2</sup> This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 19691).

### References

1. Kondo, K., Ogawa, H., Yamashita, H., *et al.* 7-Chloro-5-hydroxy-1-[2-methyl-4-(2-methylbenzoyl-amino)benzoyl]-2,3,4,5-tetrahydro-1H-1-benzazepine (OPC-41061): A potent, orally active nonpeptide arginine vasopressin V<sub>2</sub> receptor antagonist. *Bioorg. Med. Chem.* **7(8)**, 1743-1754 (1999).
2. Cadwallader, A.B., de la Torre, X., Tieri, A., *et al.* The abuse of diuretics as performance-enhancing drugs and masking agents in sport doping: Pharmacology, toxicology and analysis. *Br. J. Pharmacol.* **161(1)**, 1-16 (2010).
3. Mazzarino, M., Buccilli, V., de la Torre, X., *et al.* Characterization of the phase I and phase II metabolic profile of tolvaptan by *in vitro* studies and liquid chromatography-mass spectrometry profiling: Relevance to doping control analysis. *J. Pharm. Biomed. Anal.* **145**, 555-568 (2017).

#### 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, 07/17/2023

#### 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