

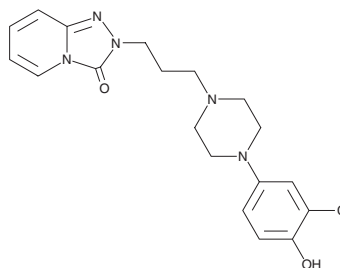
PRODUCT INFORMATION



4'-hydroxy Trazodone

Item No. 34409

CAS Registry No.: 53818-10-3
Formal Name: 2-[3-[4-(3-chloro-4-hydroxyphenyl)-1-piperazinyl]propyl]-1,2,4-triazolo[4,3-a]pyridin-3(2H)-one
MF: C₁₉H₂₂ClN₅O₂
FW: 387.9
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

4'-hydroxy Trazodone is supplied as a solid. A stock solution may be made by dissolving the 4'-hydroxy trazodone in the solvent of choice, which should be purged with an inert gas. 4'-hydroxy Trazodone is slightly soluble in chloroform and methanol (sonicated).

Description

4'-hydroxy Trazodone is a metabolite of the antidepressant and sedative trazodone.¹ It is an inhibitor of organic anion transporter 3 (OAT3; K_i = 16.9 μM) and is selective for OAT3 over OAT1 (K_i = >200 μM).²

References

1. Yamato, C., Takahashi, T., Fujita, T., *et al.* Studies on metabolism of trazodone, II. Metabolic fate after intravenous administration and effects on liver microsomal drug-metabolizing enzymes in rats. *Xenobiotica* **4**(12), 765-777 (1974).
2. Zou, L., Matsson, P., Stecula, A., *et al.* Drug metabolites potently inhibit renal organic anion transporters, OAT1 and OAT3. *J. Pharm. Sci.* **110**(1), 347-353 (2021).

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, 11/14/2022

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