

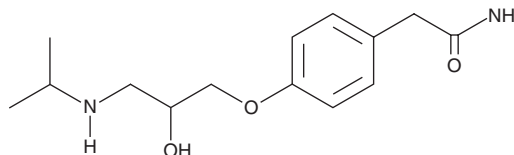
PRODUCT INFORMATION



Atenolol

Item No. 26284

CAS Registry No.: 29122-68-7
Formal Name: 4-[2-hydroxy-3-[(1-methylethyl)amino]propoxy]-benzeneacetamide
Synonyms: Duraatenol, ICI 66082
MF: C₁₄H₂₂N₂O₃
FW: 266.3
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 228, 277, 283 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

Atenolol (Item No. 26284) is an analytical reference standard categorized as a β_1 -adrenergic receptor antagonist.¹ Formulations containing atenolol have been used to enhance physical performance in athletes.² This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 17250).

References

1. Mehvar, R. and Brocks, D.R. Stereospecific pharmacokinetics and pharmacodynamics of β -adrenergic blockers in humans. *J. Pharm. Pharm. Sci.* **4**(2), 185-200 (2001).
2. Gupta, P. and Goyal, R.N. Amino functionalized graphene oxide and polymer nanocomposite based electrochemical platform for sensitive assay of anti-doping drug atenolol in biological fluids. *J. Electrochem. Soc.* **163**(13), B601-B608 (2016).

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, 07/19/2023

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