

PRODUCT INFORMATION



Dorzolamide (hydrochloride)

Item No. 26303

CAS Registry No.: 130693-82-2
Formal Name: (4S,6S)-4-(ethylamino)-5,6-dihydro-6-methyl-4H-thieno[2,3-b]thiopyran-2-sulfonamide, 7,7-dioxide, monohydrochloride

Synonyms: L-671,152, MK-507

MF: C₁₀H₁₆N₂O₄S₃ • HCl

FW: 360.9

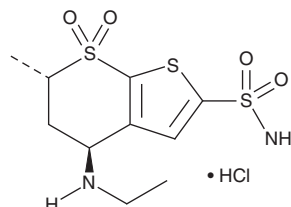
Purity: ≥98%

UV/Vis.: λ_{max}: 255 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

Dorzolamide (hydrochloride) (Item No. 26303) is an analytical reference standard categorized as a carbonic anhydrase inhibitor diuretic.¹ Diuretics have been used in sports doping to reduce weight or anabolic androgenic steroid-induced water retention, or as masking agents.² This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 14616).

References

1. Mottram, D.R. Diuretics and other masking agents. *Drugs in Sport*. Mottram, D.R., editor, 5th edition, Routledge (2011).
2. Morra, V., Davit, P., Capra, P., et al. Fast gas chromatographic/mass spectrometric determination of diuretics and masking agents in human urine: Development and validation of a productive screening protocol for antidoping analysis. *J. Chromatogr. A* **1135**(2), 219-229 (2006).

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