

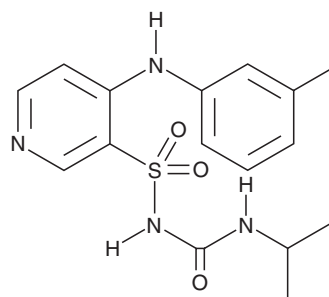
PRODUCT INFORMATION



Torsemide

Item No. 21312

CAS Registry No.: 56211-40-6
Formal Name: N-[[[(1-methylethyl)amino]carbonyl]-4-[(3-methylphenyl)amino]-3-pyridinesulfonamide
Synonyms: BM 02015, Torsemide
MF: C₁₆H₂₀N₄O₃S
FW: 348.4
Purity: ≥98%
UV/Vis.: λ_{max}: 287 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

Torsemide (Item No. 21312) is an analytical reference standard categorized as a diuretic.¹⁻³ Diuretics, including torsemide, have been abused as performance-enhancing drugs and masking agents in doping in sports.³ This product is intended for research and forensic applications.

References

1. Reyes, A.J. Diuretics in the therapy of hypertension. *J. Hum. Hypertens.* **16**, 78-83 (2002).
2. Beck, F.-X., Sone, M., Dorge, A., *et al.* Effect of loop diuretics on organic osmolytes and cell electrolytes in the renal outer medulla. *Kidney Int.* **42(4)**, 843-850 (1992).
3. 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).

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, 07/06/2023

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