

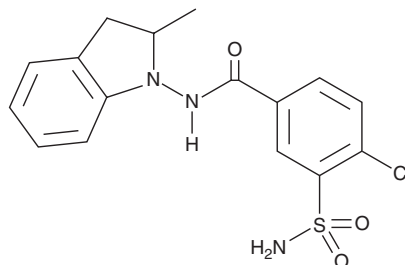
PRODUCT INFORMATION



Indapamide

Item No. 21308

CAS Registry No.: 26807-65-8
Formal Name: 3-(aminosulfonyl)-4-chloro-N-(2,3-dihydro-2-methyl-1H-indol-1-yl)-benzamide
Synonym: (±)-Indapamide
MF: C₁₆H₁₆ClN₃O₃S
FW: 365.8
Purity: ≥98%
UV/Vis.: λ_{max}: 242, 286 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

Indapamide (Item No. 21308) is an analytical reference standard that is categorized as a sulfonamide diuretic that blocks delayed-rectifier potassium currents.¹ Formulations containing indapamide are used to treat hypertension, but it is abused by athletes to reduce body weight rapidly or mask the presence of other athletic-enhancing substances.^{2,3} Indapamide can be detected in urine.⁴ This product is intended for research and forensic applications.

References

1. Turgeon, J., Daleau, P., Bennett, P.B., *et al.* Block of IKS, the slow component of the delayed rectifier K⁺ current, by the diuretic agent indapamide in guinea pig myocytes. *Circ. Res.* **75(5)**, 879-886 (2017).
2. Robinson, D.M. and Wellington, K. Indapamide sustained release: A review of its use in the treatment of hypertension. *Drugs* **66(2)**, 257-271 (2006).
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).
4. Ramkumar, A., Ponnusamy, V.K., and Jen, J.-F. Rapid determination of indapamide in human urine using novel low-density solvent based ultrasound assisted emulsification microextraction coupled with high performance liquid chromatography-variable wavelength detection. *Anal. Methods* **5**, 2572-2577 (2013).

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, 06/28/2023

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