

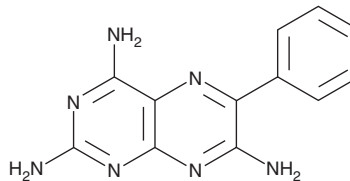
PRODUCT INFORMATION



Triamterene

Item No. 22486

CAS Registry No.: 396-01-0
Formal Name: 6-phenyl-2,4,7-pteridinetriamine
Synonyms: NSC 77625, SKF 8542,
2,4,7-Triamino-6-phenylpteridine
MF: C₁₂H₁₁N₇
FW: 253.3
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 266, 366 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

Triamterene (Item No. 22486) is an analytical reference standard that is categorized as a pteridine. Triamterene is an inhibitor of the epithelial sodium channel (ENaC) that increases renal sodium excretion and decreases potassium excretion. It inhibits rat ENaC with an IC₅₀ value of 4.5 μM.¹ Diuretics, including triamterene, have been abused as performance-enhancing drugs and masking agents in sports doping.² This product is intended for research and forensic applications.

References

1. Kellenberger, S., Gautschi, I., and Schlid, L. Mutations in the epithelial Na⁺ channel ENaC outer pore disrupt amiloride block by increasing its dissociation rate. *Mol. Pharmacol.* **64(4)**, 848-856 (2003).
2. 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, 06/13/2023

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