

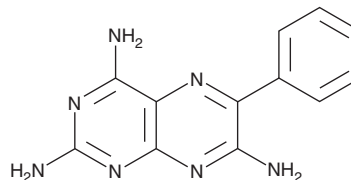
PRODUCT INFORMATION



Triamterene

Item No. 21242

CAS Registry No.: 396-01-0
Formal Name: 6-phenyl-2,4,7-pteridinetriamine
Synonyms: NSC 77625, SKF 8542
MF: C₁₂H₁₁N₇
FW: 253.3
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 271, 367 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Triamterene is supplied as a crystalline solid. A stock solution may be made by dissolving the triamterene in the solvent of choice, which should be purged with an inert gas. Triamterene is soluble in the organic solvent ethanol. The solubility of triamterene in ethanol is approximately 1 mg/ml.

Triamterene is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, triamterene should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Triamterene has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Triamterene is an inhibitor of the epithelial sodium channel (ENaC; IC₅₀ = 4.5 μM for the rat channel).¹ *In vivo*, triamterene (0.5-32 mg/animal) enhances sodium secretion and decreases potassium secretion in adrenalectomized rats.² Formulations containing triamterene have been used in the treatment of edema. This product is also available as an analytical reference standard (Item No. 22486).

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. Baba, W.I., Tudhope, G.R., and Wilson, G.M. Triamterene, a new diuretic drug. I. Studies in normal men and in adrenalectomized rats. *Br. Med. J.* **2**(5307), 756-760 (1962).

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, 10/18/2022

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