

PRODUCT INFORMATION



Azaprocin (hydrochloride)

Item No. 42268

Formal Name: 8-(1-oxopropyl)-3-[(2E)-3-phenyl-2-propenyl]-3,8-diazabicyclo[3.2.1]octane, monohydrochloride

MF: C₁₈H₂₄N₂O • HCl

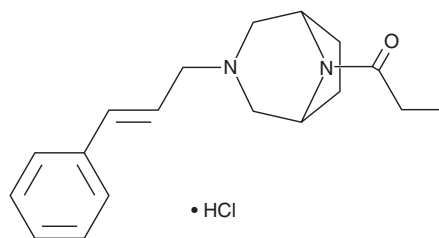
FW: 320.9

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Description

Azaprocin (hydrochloride) (Item No. 42268) is an analytical reference standard categorized as an opioid.^{1,2} Azaprocin has analgesic activity in rodents.¹ This product is intended for research and forensic applications.

References

1. Cignarella, G., Occelli, E., Maffii, G., *et al.* Bicyclic homologs of piperazine. VII. Synthesis of pharmacologically active 8-methyl-3,8-diazabicyclo(3.2.1)octanes. III. *J. Med. Chem.* **6**, 29-36 (1963).
2. Cignarella, G., Barlocco, D., Vianello, P., *et al.* Benzocondensed derivatives as rigid analogues of the μ -opioid agonist 3(8)-cinnamyl-8(3)-propionyl-3,8-diazabicyclo[3.2.1]octanes: synthesis, modeling, and affinity. *ChemMedChem* **53**, 667-674 (1998).

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, 09/16/2025

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