

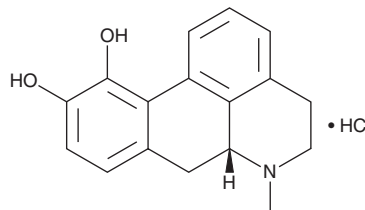
PRODUCT INFORMATION



(-)-Apomorphine (hydrochloride)

Item No. 34222

CAS Registry No.: 314-19-2
Formal Name: 5,6,6aR,7-tetrahydro-6-methyl-4H-dibenzo[de,g]quinoline-10,11-diol, monohydrochloride
Synonym: NSC 11442
MF: C₁₇H₁₇NO₂ • HCl
FW: 303.8
Purity: ≥98%
UV/Vis.: λ_{max}: 274 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

(-)-Apomorphine (hydrochloride) (Item No. 34222) is an analytical reference standard categorized as a dopamine receptor agonist.¹ This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 16094).

Reference

1. Millan, M.J., Maiorini, L., Cussac, D., *et al.* Differential actions of antiparkinson agents at multiple classes of monoaminergic receptor. I. A multivariate analysis of the binding profiles of 14 drugs at 21 native and cloned human receptor subtypes. *J. Pharmacol. Exp. Ther.* **303**(2), 791-804 (2002).

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