

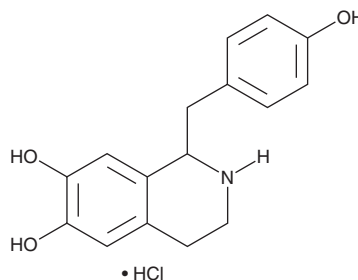
PRODUCT INFORMATION



(±)-Higenamine (hydrochloride)

Item No. 24792

CAS Registry No.: 11041-94-4
Formal Name: 1,2,3,4-tetrahydro-1-[(4-hydroxyphenyl)methyl]-6,7-isoquinolinediol, monohydrochloride
Synonyms: (±)-Demethylcoclaurine, Norcoclaurine
MF: C₁₆H₁₇NO₃ • HCl
FW: 307.8
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

(±)-Higenamine (hydrochloride) (Item No. 24792) is an analytical reference standard categorized as a β -adrenergic receptor agonist.¹ (±)-Higenamine has been found in performance-enhancing drug (PED) samples seized by law enforcement.² This product is intended for research and forensic applications.

References

1. Chen, Y., Guo, B., Zhang, H., *et al.* Higenamine, a dual agonist for β_1 - and β_2 -adrenergic receptors identified by screening a traditional Chinese medicine library. *Planta Med.* 85(9-10), 738-744 (2019).
2. Grucza, K., Kowalczyk, K., Wicka, M., *et al.* The use of a valid and straightforward method for the identification of higenamine in dietary supplements in view of anti-doping rule violation cases. *Drug Test Anal.* 11(6), 912-917 (2019).

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, 08/19/2024

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