

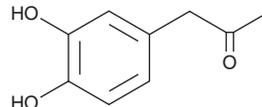
PRODUCT INFORMATION



3',4'-Dihydroxyphenylacetone

Item No. 19522

CAS Registry No.: 2503-44-8
Formal Name: 1-(3,4-dihydroxyphenyl)-2-propanone
MF: C₉H₁₀O₃
FW: 166.2
Purity: ≥98%
UV/Vis.: λ_{max}: 284 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.

Description

3',4'-Dihydroxyphenylacetone (Item No. 19522) is an analytical reference standard. It is a minor metabolite of 3,4-MDEA (Item Nos. 14085 | 15689), MDMA, and α-methyl dopa produced by oxidative deamination.¹⁻³ This product is intended for forensic and research applications.

References

1. Lim, H.K. and Foltz, R.L. *In vivo* and *in vitro* metabolism of 3,4-(methylenedioxy)methamphetamine in the rat: Identification of metabolites using an ion trap detector. *Chem. Res. Toxicol.* **1(6)**, 370-378 (1988).
2. Ensslin, H.K., Maurer, H.H., Gouzoulis, E., *et al.* Metabolism of racemic 3,4-methylenedioxyethylamphetamine in humans. Isolation, identification, quantification, and synthesis of urinary metabolites. *Drug Metab. Dispos.* **24(8)**, 813-820 (1996).
3. Bertoldi, M., Dominici, P., Moore, P.S., *et al.* Reaction of dopa decarboxylase with α-methyl dopa leads to an oxidative deamination producing 3,4-dihydroxyphenylacetone, an active site directed affinity label. *Biochem.* **37(18)**, 6552-6561 (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, 11/23/2022

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