

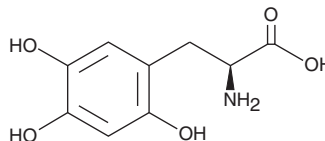
PRODUCT INFORMATION



6-hydroxy-L-DOPA

Item No. 45703

CAS Registry No.: 27244-64-0
Formal Name: 2,5-dihydroxy-L-tyrosine
Synonyms: L-6-hydroxy DOPA,
Levodopa Related Compound A
MF: $C_9H_{11}NO_5$
FW: 213.2
Purity: $\geq 95\%$
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.

Laboratory Procedures

6-hydroxy-L-DOPA is supplied as a solid. Aqueous solutions of 6-hydroxy-L-DOPA can be prepared by directly dissolving the solid in aqueous buffers. 6-hydroxy-L-DOPA is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

6-hydroxy-L-DOPA is a derivative of L-DOPA (Item No. 13248).¹ It has been used as an electron-donating unit in luminescent silver (Ag) nanocluster sensors for the detection of nitro-substituted explosives.

Reference

1. Enkin, N., Sharon, E., Golub, E., *et al.* Ag nanocluster/DNA hybrids: Functional modules for the detection of nitroaromatic and RDX explosives. *Nano Lett.* **14**(8), 4918-4922 (2014).

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, 06/22/2026

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