

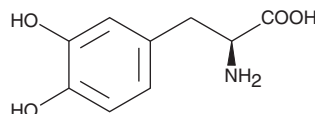
# PRODUCT INFORMATION



## L-DOPA

Item No. 13248

**CAS Registry No.:** 59-92-7  
**Formal Name:** 3-hydroxy-L-tyrosine  
**Synonyms:** 3,4-Dihydroxyphenylalanine, Levodopa  
**MF:** C<sub>9</sub>H<sub>11</sub>NO<sub>4</sub>  
**FW:** 197.2  
**Purity:** ≥95%  
**Supplied as:** A crystalline solid  
**Storage:** Room temperature  
**Stability:** ≥4 years



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

### Laboratory Procedures

L-DOPA is supplied as a crystalline solid. L-DOPA is sparingly soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

L-DOPA is a metabolic precursor of dopamine that is capable of crossing the blood brain barrier.<sup>1</sup> It is produced from L-tyrosine by tyrosine hydroxylase and metabolized by catechol-O-methyl transferase (COMT). In the brain L-DOPA is converted to dopamine. Formulations containing L-DOPA have been used to increase dopamine concentrations in the brain as a treatment for Parkinson's disease and stroke recovery.

### Reference

1. Berends, H.I., Nijlant, J.M.M., Movig, K.L.L., *et al.* The clinical use of drugs influencing neurotransmitters in the brain to promote motor recovery after stroke; a Cochrane systematic review. *Eur J Phys Rehabil Med* 45(4), 621-630 (2009).

#### 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, 12/19/2022

#### 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