

# PRODUCT INFORMATION

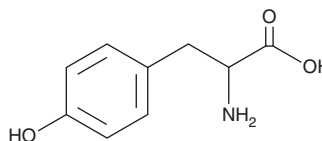


## DL-Tyrosine

Item No. 36460

CAS Registry No.: 556-03-6  
Synonyms: DL-4-Hydroxyphenylalanine,  
DL-*p*-Hydroxyphenylalanine,  
DL-*para*-Hydroxyphenylalanine,  
NSC 205006, (±)-Tyrosine

MF: C<sub>9</sub>H<sub>11</sub>NO<sub>3</sub>  
FW: 181.2  
Purity: ≥98%  
UV/Vis.: λ<sub>max</sub>: 226, 282 nm  
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

DL-Tyrosine is supplied as a solid. A stock solution may be made by dissolving the DL-tyrosine in the solvent of choice, which should be purged with an inert gas. DL-Tyrosine is slightly soluble in DMSO.

DL-Tyrosine is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

DL-Tyrosine is a racemic mixture of the atypical amino acid D-tyrosine and the conditionally essential amino acid L-tyrosine (Item No. 36333).<sup>1,2</sup>

### References

1. Grishin, D.V., Zhdanov, D.D., Pokrovskaya, M.V., *et al.* D-amino acids in nature, agriculture and biomedicine. *All Life* **13(1)**, 11-22 (2020).
2. Kopple, J.D. Phenylalanine and tyrosine metabolism in chronic kidney failure. *J. Nutr.* **137(6 Suppl 1)**, 1586S-1590S (2007).

#### 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, 04/20/2023

#### 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