

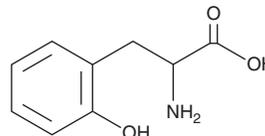
PRODUCT INFORMATION



DL-*o*-Tyrosine

Item No. 33570

CAS Registry No.: 2370-61-8
Formal Name: 2-hydroxy-phenylalanine
Synonyms: DL-2-Hydroxyphenylalanine, NSC 72345
MF: C₉H₁₁NO₃
FW: 181.2
Purity: ≥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

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

Description

DL-*o*-Tyrosine is a hydroxylation product of phenylalanine.¹ It is formed from phenylalanine by hydroxyl radicals generated during radiolysis and can be further hydroxylated to form 2,3-dihydroxyphenylalanine (2,3-DOPA). DL-*o*-Tyrosine has been used as a marker of irradiation processing in poultry meat.²

References

1. Żegota, H., Kołodziejczyk, K., Król, M., *et al.* *o*-Tyrosine hydroxylation by OH• radicals. 2,3-DOPA and 2,5-DOPA formation in γ -irradiated aqueous solution. *Radiat. Phys. Chem.* **72(1)**, 25-33 (2005).
2. Chuaqui-Offermans, N. and McDougall, T. Background levels and radiation dose yield of *o*-tyrosine in chicken meat. *J. Food Prot.* **54(12)**, 935-938 (1991).

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, 11/15/2022

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