

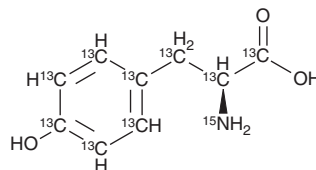
PRODUCT INFORMATION



L-Tyrosine-¹³C₉,¹⁵N

Item No. 31256

CAS Registry No.: 202407-26-9
Formal Name: L-tyrosine-1,2,3,4,5,6-¹³C₉-¹⁵N
Synonyms: L-4-Hydroxyphenylalanine,
p-Hydroxyphenylalanine,
(-)-Tyrosine, p-Tyrosine
MF: [¹³C]₉H₁₁[¹⁵N]O₃
FW: 191.1
Purity: ≥98%
UV/Vis.: λ_{max}: 223 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

L-Tyrosine-¹³C₉,¹⁵N is supplied as a solid. A stock solution may be made by dissolving the L-tyrosine-¹³C₉,¹⁵N in the solvent of choice, which should be purged with an inert gas. L-Tyrosine-¹³C₉,¹⁵N is slightly soluble in methanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of L-tyrosine-¹³C₉,¹⁵N can be prepared by directly dissolving the L-tyrosine-¹³C₉,¹⁵N in aqueous buffers. L-Tyrosine-¹³C₉,¹⁵N is slightly soluble in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

L-Tyrosine-¹³C₉,¹⁵N is intended for use as an internal standard for the quantification of L-tyrosine by GC- or LC-MS. L-Tyrosine is a conditionally essential amino acid.¹ It is produced by hydroxylation of phenylalanine by phenylalanine hydroxylase but can also be obtained from dietary sources or degradation of endogenous proteins, resulting in L-tyrosine release.^{1,2} L-Tyrosine is a precursor in the biosynthesis of catecholamine neurotransmitters, melanins, and thyroid hormones.³ Plasma, skeletal muscle, and erythrocyte levels of L-tyrosine are decreased in patients with chronic kidney disease.¹

References

1. Kopple, J.D. Phenylalanine and tyrosine metabolism in chronic kidney failure. *J. Nutr.* **137(6 Suppl 1)**, 1586S-1590S (2007).
2. Webster, D. and Wildgoose, J. Tyrosine supplementation for phenylketonuria. *Cochrane DB Syst. Rev.* **6**, CD001507 (2013).
3. Slominski, A. and Paus, R. Towards defining receptors for L-tyrosine and L-DOPA. *Mol. Cell Endocrinol.* **99(2)**, C7-C11 (1994).

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, 04/19/2023

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