

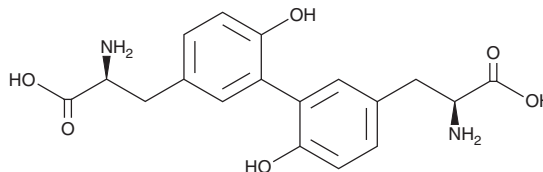
# PRODUCT INFORMATION



## L,L-Dityrosine

Item No. 35746

**CAS Registry No.:** 63442-81-9  
**Formal Name:**  $\alpha^3S, \alpha^3'S$ -diamino-6,6'-dihydroxy-[1,1'-biphenyl]-3,3'-dipropanoic acid  
**MF:**  $C_{18}H_{20}N_2O_6$   
**FW:** 360.4  
**Purity:**  $\geq 90\%$   
**Supplied as:** A solution in ethanol  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 2$  years



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

### Description

L,L-Dityrosine is an amino acid conjugate containing two L-tyrosine molecules. L-Tyrosine residues in fibrinogen proteins can form L,L-dityrosine cross-links between fibrinogen proteins in the presence of hematin and plasma-generated hydrogen peroxide under non-thermal plasma exposure.<sup>1</sup> L-Tyrosine residues in human, but not mouse, amylin peptides can form L,L-dityrosine cross-links between amylin peptides in the presence of copper (II) ions and hydrogen peroxide.<sup>2</sup>

### References

1. Ke, Z. and Huang, Q. Haem-assisted dityrosine-cross-linking of fibrinogen under non-thermal plasma exposure: One important mechanism of facilitated blood coagulation. *Sci. Rep.* **6**, 26982 (2016).
2. Dong, X., Svantesson, T., Sholts, S.B., *et al.* Copper ions induce dityrosine-linked dimers in human but not in murine islet amyloid polypeptide (IAPP/amylin). *Biochem. Biophys. Res. Commun.* **510(4)**, 520-524 (2019).

#### 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, 11/09/2022

#### 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