

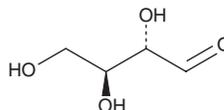
PRODUCT INFORMATION



L-(+)-Erythrose

Item No. 20381

CAS Registry No.: 533-49-3
Formal Name: (2S,3S)-2,3,4-trihydroxy-butanal
Synonyms: (+)-Erythrose, L-Erythrose
MF: C₄H₈O₄
FW: 120.1
Purity: ≥95%
Supplied as: A 0.108 M aqueous solution
Storage: -20°C
Stability: ≥2 years



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

Description

L-(+)-Erythrose is an aldotetrose carbohydrate that has been used in glycation studies and to characterize erythrose reductase activity.^{1,2}

References

1. Li, Y., Drew, M.G.B., Welchman, E.V., *et al.* Syntheses of ethyl 3-deoxy-3,3-difluoro-D-arabino-heptulosonate and analogues. *Tetrahedron* **60(31)**, 6523-6531 (2004).
2. Del-Corso, A., Balestri, F., Di Bugno, E., *et al.* A new approach to control the enigmatic activity of aldose reductase. *PLoS One* **8(9)**, (2013).

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, 12/13/2022

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