

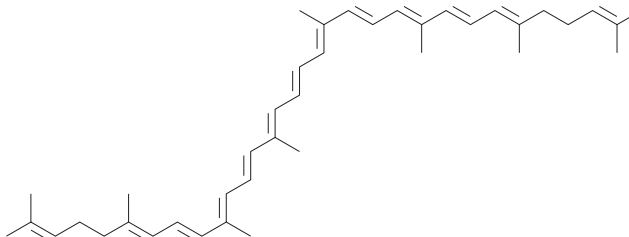
PRODUCT INFORMATION



Lycopene

Item No. 70945

CAS Registry No.: 502-65-8
Formal Name: ψ,ψ -carotene
Synonym: NSC 407322
MF: $C_{40}H_{56}$
FW: 536.9
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: -80°C
Stability: ≥ 2 years
Special Conditions: Light/temperature sensitive, not stable in solution.



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

Laboratory Procedures

Lycopene is supplied as a crystalline solid. A stock solution may be made by dissolving the lycopene in the solvent of choice. Lycopene is soluble in chloroform at a concentration of approximately 3 mg/ml.

Lycopene is sparingly 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

Lycopene is a red-colored carotenoid found in tomatoes and other red fruits and vegetables.¹ Carotenoids, including lycopene, are powerful antioxidants that efficiently quench singlet oxygen.^{1,2} Presumably through this action, carotenoids may protect against cancers, cardiovascular stress, and other diseases.³⁻⁶

References

1. Stahl, W., Heinrich, U., Aust, O., *et al.* Lycopene-rich products and dietary photoprotection. *Photochem. Photobiol. Sci.* **5**(2), 238-242 (2006).
2. Cantrell, A., McGarvey, D.J., Truscott, T.G., *et al.* Singlet oxygen quenching by dietary carotenoids in a model membrane environment. *Arch. Biochem. Biophys.* **412**(1), 47-54 (2003).
3. Vance, T.M., Su, J., Fontham, E.T., *et al.* Dietary antioxidants and prostate cancer: A review. *Nutr. Cancer* **65**(6), 793-801 (2013).
4. Sato, R., Helzlsouer, K.J., Alberg, A.J., *et al.* Prospective study of carotenoids, tocopherols, and retinoid concentrations and the risk of breast cancer. *Cancer Epidemiol. Biomarkers Prev.* **11**(5), 451-457 (2002).
5. Giordano, P., Scicchitano, P., Locorotondo, M., *et al.* Carotenoids and cardiovascular risk. *Curr. Pharm. Des.* **18**(34), 5577-5589 (2012).
6. Aune, D., Chan, D.S.M., Vieira, A.R., *et al.* Dietary compared with blood concentrations of carotenoids and breast cancer risk: A systematic review and meta-analysis of prospective studies. *Am. J. Clin. Nutr.* **96**(2), 356-373 (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, 12/21/2023

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