

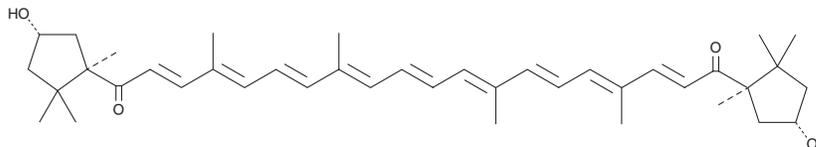
PRODUCT INFORMATION



Capsorubin

Item No. 35115

CAS Registry No.: 470-38-2
Formal Name: (3S,3'S,5R,5'R)-3,3'-dihydroxy-
κ,κ-carotene-6,6'-dione
MF: C₄₀H₅₆O₄
FW: 600.9
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Capsorubin is supplied as a solid. A stock solution may be made by dissolving the capsorubin in the solvent of choice, which should be purged with an inert gas. Capsorubin is soluble in DMSO and dimethyl formamide.

Description

Capsorubin is a carotenoid that has been found in *C. annuum* and has diverse biological activities.¹⁻³ It inhibits lipid peroxidation induced by 2,2'-azobis(2,4-dimethyl valeronitrile) (AMVN) *in vitro* when used at a concentration of 167 μM.¹ Capsorubin (1 μM) decreases UVB-induced formation of DNA strand breaks and apoptosis in isolated human dermal fibroblasts.² It also inhibits Epstein-Barr virus early antigen (EBV-EA) activation induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014), a marker of tumor formation, in Raji cells.

References

1. Maoka, T., Goto, Y., Isobe, K., *et al.* Antioxidative activity of capsorubin and related compounds from paprika (*Capsicum annuum*). *J. Oleo Sci.* **50(8)**, 663-665 (2001).
2. Fernández-García, E., Carvajal-Lérida, I., and Pérez-Gálvez, A. Carotenoids exclusively synthesized in red pepper (capsanthin and capsorubin) protect human dermal fibroblasts against UVB induced DNA damage. *Photochem. Photobiol. Sci.* **15(9)**, 1204-1211 (2016).
3. Maoka, T., Mochida, K., Kozuka, M., *et al.* Cancer chemopreventive activity of carotenoids in the fruits of red paprika *Capsicum annuum* L. *Cancer Lett.* **172(2)**, 103-109 (2001).

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, 11/21/2022

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