

PRODUCT INFORMATION

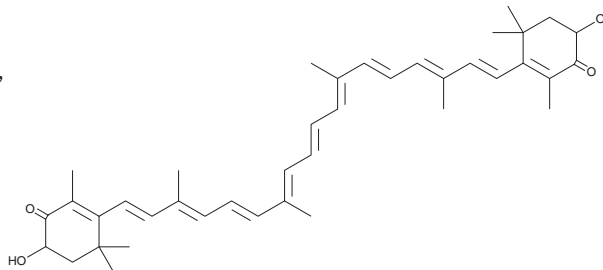


rac-Astaxanthin

Item No. 70685

CAS Registry No.: 7542-45-2
Formal Name: 3,3'-dihydroxy- β,β -carotene-4,4'-dione
Synonyms: AstaREAL, AstaXin, BioAstin, Carophyll Pink, Lucantin Pink, NatuRose, Ovoester

MF: C₄₀H₅₂O₄
FW: 596.9
Purity: \geq 97%
UV/Vis.: λ_{max} : 473 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

rac-Astaxanthin is supplied as a crystalline solid. A stock solution may be made by dissolving the *rac*-astaxanthin in the solvent of choice, which should be purged with an inert gas. *rac*-Astaxanthin is soluble in organic solvents such as ethanol, chloroform, DMSO, and dimethyl formamide. The solubility of *rac*-astaxanthin in these solvents is approximately 0.09, 6, 0.13, and 0.5 mg/ml, respectively. *rac*-Astaxanthin is extremely lipophilic and crashes out of solution (of organic solvents), even at the smallest addition of water.

Description

rac-Astaxanthin is a carotenoid pigment found primarily in marine animals including shrimp and salmon.¹ It is a potent lipid-soluble antioxidant.

Reference

- Schroeder, W.A. and Johnson, E.A. Singlet oxygen and peroxy radicals regulate carotenoid biosynthesis in *Phaffia rhodozyma*. *J. Biol. Chem.* **270**(31), 18374-18379 (1995).

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, 04/06/2021

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