

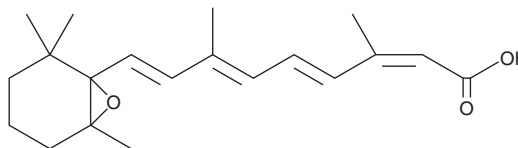
PRODUCT INFORMATION



5,6-epoxy-13-cis Retinoic Acid

Item No. 23091

CAS Registry No.: 81444-57-7
Formal Name: 5,6-epoxy-5,6-dihydro-13-cis-retinoic acid
Synonym: 5,6-epoxy-13-cis RA
MF: C₂₀H₂₈O₃
FW: 316.4
Purity: ≥95%
Supplied as: A solid
Storage: -80°C
Stability: ≥2 years



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

Laboratory Procedures

5,6-epoxy-13-cis Retinoic acid is supplied as a solid. A stock solution may be made by dissolving the 5,6-epoxy-13-cis retinoic acid in the solvent of choice, which should be purged with an inert gas. 5,6-epoxy-13-cis Retinoic acid is slightly soluble in methanol (sonicated).

Description

5,6-epoxy-13-cis Retinoic acid is a metabolite of 13-cis retinoic acid.¹ It is formed when 13-cis retinoic acid undergoes cooxidation by prostaglandin H (PGH) synthase in the presence of hydroperoxides or peroxy radicals. 5,6-epoxy-13-cis Retinoic acid is also a potential impurity found in commercial preparations of 13-cis retinoic acid.²

References

1. Samokyszyn, V.M. and Marnett L.J. Hydroperoxide-dependent cooxidation of 13-cis-retinoic acid by prostaglandin H synthase. *J. Biol. Chem.* **262(29)**, 14119-14133 (1987).
2. Madan, H.K., Venkateshwaran, R., Madan, S., *et al.* Low dose oral pharmaceutical composition of isotretinoin. *US9750711B2* (2017).

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, 02/19/2021

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