

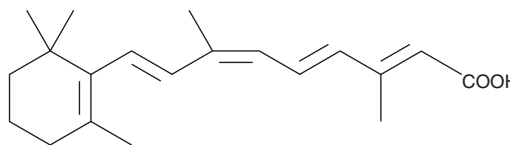
PRODUCT INFORMATION



9-*cis*-Retinoic Acid

Item No. 14587

CAS Registry No.: 5300-03-8
Formal Name: 9-*cis*-retinoic acid
Synonyms: Alitretinoin, NSC 659772, Panretin, 9-*cis*-RA
MF: C₂₀H₂₈O₂
FW: 300.4
Purity: ≥95%
UV/Vis.: λ_{max}: 42, 348, 357 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

9-*cis*-Retinoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 9-*cis*-retinoic acid in the solvent of choice. 9-*cis*-Retinoic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of 9-*cis*-retinoic acid in ethanol is approximately 0.5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

9-*cis*-Retinoic acid 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

9-*cis*-Retinoic acid is a natural metabolite of vitamin A, derived from the intermediate all-*trans* retinoic acid (Item No. 11017).¹ It potently activates all isoforms of retinoic acid receptor (RAR) (K_i = 0.5-27 nM) as well as retinoid X receptor (RXR) isoforms (K_i = 3.8-12 nM).^{2,3} RAR heterodimerizes with RXR, while RXR can homodimerize as well as heterodimerize with numerous partners in addition to RAR, thus allowing 9-*cis*-retinoic acid to evoke a wide range of effects.^{1,4}

References

1. Kane, M.A. Analysis, occurrence, and function of 9-*cis*-retinoic acid. *Biochim. Biophys. Acta* **1821**(1), 10-20 (2012).
2. Wong, M.F., Repa, J.J., Clagett-Dame, M., et al. Synthesis and receptor binding affinity of conformationally restricted retinoic acid analogues. *Bioorg. Med. Chem. Lett.* **7**(17), 2313-2318 (1997).
3. Umemiya, H., Fukasawa, H., Ebisawa, M., et al. Regulation of retinoid actions by diazepinylbenzoic acids. Retinoid synergists which activate the RXR-RAR heterodimers. *J. Med. Chem.* **40**(26), 4222-4234 (1997).
4. Dawson, M.I. and Xia, Z. The retinoid X receptors and their ligands. *Biochim Biophys. Acta.* **1821**(1), 21-56 (2012).

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, 09/19/2017

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