

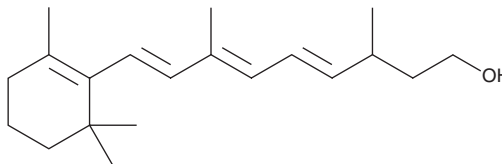
PRODUCT INFORMATION



all-*trans*-13,14-Dihydroretinol

Item No. 34873

CAS Registry No.: 115797-14-3
Formal Name: 13,14-dihydro-retinol
MF: $C_{20}H_{32}O$
FW: 288.5
Purity: $\geq 80\%$
Supplied as: A 5 mg/ml solution in ethyl acetate
Storage: -80°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

all-*trans*-13,14-Dihydroretinol is supplied as a solution in ethyl acetate. To change the solvent, simply evaporate the ethyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. all-*trans*-13,14-Dihydroretinol is slightly soluble in chloroform and methanol.

Description

all-*trans*-13,14-Dihydroretinol is a metabolite of all-*trans* retinoic acid (Item No. 11017).¹ all-*trans*-13,14-Dihydroretinol is formed from all-*trans* retinoic acid by retinol saturase.

Reference

1. Weber, P., Flores, R.E., Kiefer, M.F., *et al.* Retinol saturase: More than the name suggests. *Trends Pharmacol. Sci.* **41**(6), 418-427 (2020).

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, 02/17/2026

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