

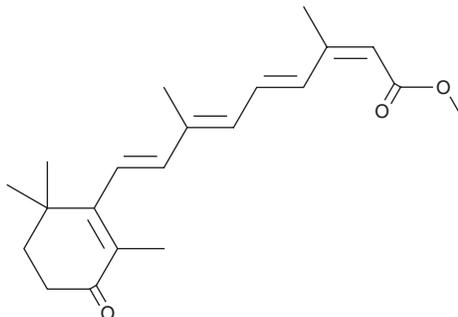
# PRODUCT INFORMATION



## Methyl 13-cis-4-oxoretinoate

Item No. 38530

**CAS Registry No.:** 71748-57-7  
**Formal Name:** 13-cis-4-oxo-retinoic acid, methyl ester  
**Synonyms:** Methyl 13-cis-4-oxo RA,  
Methyl 13-cis-4-oxo Retinoic Acid  
**MF:** C<sub>21</sub>H<sub>28</sub>O<sub>3</sub>  
**FW:** 328.5  
**Purity:** ≥85%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Methyl 13-cis-4-oxoretinoate is supplied as a solid. A stock solution may be made by dissolving the methyl 13-cis-4-oxoretinoate in the solvent of choice, which should be purged with an inert gas. Methyl 13-cis-4-oxoretinoate is sparingly soluble (1-10 mg/ml) in chloroform and acetonitrile.

### Description

Methyl 13-cis-4-oxoretinoate is a degradation product of 13-cis retinoic acid (isotretinoin; Item No. 21648).<sup>1</sup> It is formed from 13-cis retinoic acid under acidic conditions.

### Reference

1. Tanaka, H., Kagechika, H., and Shudo, K. Specific oxidation of retinoic acid to 4-oxo-retinoic acid in diluted acid solutions. *Chem. Pharm. Bull.* **43**(2), 356-358 (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, 09/04/2025

#### 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