

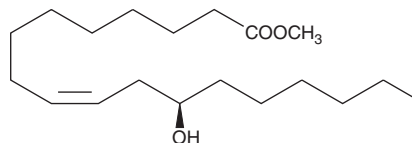
# PRODUCT INFORMATION



## Ricinoleic Acid methyl ester

Item No. 10009729

**CAS Registry No.:** 141-24-2  
**Formal Name:** 12R-hydroxy-9Z-octadecenoic acid, methyl ester  
**Synonyms:** NSC 1254, SFE 19:1;O  
**MF:** C<sub>19</sub>H<sub>36</sub>O<sub>3</sub>  
**FW:** 312.5  
**Purity:** ≥98%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

Ricinoleic acid methyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of ricinoleic acid methyl ester in these solvents is approximately 50 mg/ml.

Ricinoleic acid methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of ricinoleic acid methyl ester should be diluted with the aqueous buffer of choice. Ricinoleic acid methyl ester has a solubility of 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Ricinoleic acid is a naturally occurring 12-hydroxy fatty acid. It constitutes about 90% of the fatty acids in castor oil. Ricinoleic acid can serve as a substrate for the synthesis of conjugated linoleic acids. It is considered safe and non-toxic as a food constituent in humans, up to a daily intake of 2.5 grams per day. In high doses, castor oil has a laxative effect. Ricinoleic acid methyl ester is a neutral, more lipophilic form of the free acid that can be used as an analytical standard for ricinoleic acid.

#### 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, 03/27/2024

#### 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