

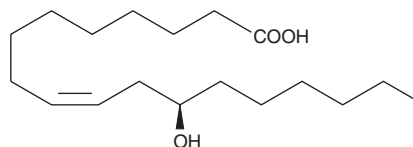
PRODUCT INFORMATION



Ricinoleic Acid

Item No. 10009369

CAS Registry No.: 141-22-0
Formal Name: 12R-hydroxy-9Z-octadecenoic acid
Synonyms: (R)-12-Hydroxyoleic Acid,
NSC 281242
MF: C₁₈H₃₄O₃
FW: 298.5
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 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of ricinoleic acid in these solvents is approximately 20 mg/ml.

Ricinoleic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of ricinoleic acid should be diluted with the aqueous buffer of choice. Ricinoleic acid 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.

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, 11/07/2022

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