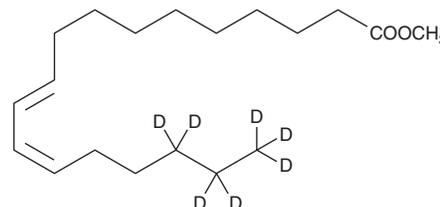


# PRODUCT INFORMATION



## 10(E),12(Z)-Conjugated Linoleic Acid-d<sub>7</sub> methyl ester Item No. 26483

<b>Formal Name:</b>	10E,12Z-octadecadienoic acid-16,16,17,17,18,18,18-d <sub>7</sub> , methyl ester
<b>Synonyms:</b>	10E,12Z-CLA-d <sub>7</sub> methyl ester, Methyl 10(E),12(Z)-Octadecadienoate-d <sub>7</sub> methyl ester, SFE 19:2-d <sub>7</sub>
<b>MF:</b>	C <sub>19</sub> H <sub>27</sub> D <sub>7</sub> O <sub>2</sub>
<b>FW:</b>	301.5
<b>Chemical Purity:</b>	≥95% (10(E),12(Z)-Conjugated Linoleic Acid methyl ester)
<b>Deuterium Incorporation:</b>	≥99% deuterated forms (d <sub>1</sub> -d <sub>7</sub> ); ≤1% d <sub>0</sub>
<b>UV/Vis.:</b>	λ <sub>max</sub> : 233 nm
<b>Supplied as:</b>	A 1 mg/ml solution in ethanol
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥2 years



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

### Laboratory Procedures

10(E),12(Z)-Conjugated linoleic acid-d<sub>7</sub> methyl ester is intended for use as an internal standard for the quantification of 10(E),12(Z)-conjugated linoleic acid methyl ester (Item No. 24584) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

10(E),12(Z)-Conjugated linoleic acid-d<sub>7</sub> 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 10(E),12(Z)-conjugated linoleic acid-d<sub>7</sub> methyl ester in these solvents is approximately 100 mg/ml.

### Description

10(E),12(Z)-Conjugated linoleic acid methyl ester is a linoleic acid methyl ester containing conjugated double bonds at the 10 and 12 positions.

**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, 12/19/2025

### 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