

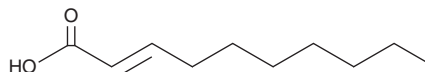
# PRODUCT INFORMATION



## *trans*-2-Decenoic Acid

Item No. 17888

**CAS Registry No.:** 334-49-6  
**Formal Name:** 2E-decenoic acid  
**Synonym:** *trans*-2-Decylenic Acid, FA 10:1  
**MF:** C<sub>10</sub>H<sub>18</sub>O<sub>2</sub>  
**FW:** 170.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 209 nm  
**Supplied as:** A neat oil  
**Storage:** -20°C  
**Stability:** ≥4 year



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

### Laboratory Procedures

*trans*-2-Decenoic acid is supplied as a neat oil. A stock solution may be made by dissolving the *trans*-2-decenoic acid in the solvent of choice, which should be purged with an inert gas. *trans*-2-Decenoic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of *trans*-2-decenoic acid ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

### Description

*trans*-2-Decenoic acid is an unsaturated fatty acid found in royal jelly produced from the hypopharyngeal and mandibular gland secretions of honeybees. It is isolated in conjunction with 10-HDA (Item No. 10976) and both demonstrate weak estrogenic activity, inhibiting binding of 17β-estradiol (Item No. ISO60155) to estrogen receptor β.<sup>1</sup>

### Reference

1. Suzuki, K.-M., Isohama, Y., Maruyama, H., *et al.* Estrogenic activities of fatty acids and sterol isolated from royal jelly. *Evid. Based Complement. Alternat. Med.* **5(3)**, 295-302 (2008).

#### 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, 02/08/2024

#### 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