

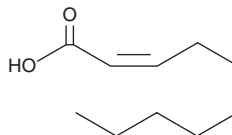
# PRODUCT INFORMATION



## *cis*-2-Decenoic Acid

Item No. 11966

CAS Registry No.: 15790-91-7  
Formal Name: 2*Z*-decenoic acid  
Synonym: FA 10:1  
MF: C<sub>10</sub>H<sub>18</sub>O<sub>2</sub>  
FW: 170.3  
Purity: ≥95%  
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

*cis*-2-Decenoic 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 ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of *cis*-2-decenoic acid in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

### Description

*cis*-2-Decenoic acid is an unsaturated short chain fatty acid that is secreted by *P. aeruginosa* and induces a dispersion response in biofilms formed by gram-negative and gram-positive bacteria, as well as by the yeast *C. albicans*.<sup>1</sup> It effectively promotes dispersion of *P. aeruginosa* biofilms over a concentration range of 1.0 to 10 nM.<sup>1</sup>

### Reference

1. Davies, D.G., and Marques, C.N.H. A fatty acid messenger is responsible for inducing dispersion in microbial biofilms. *J. Bacteriol.* **191**(5), 1393-1403 (2009).

#### 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, 02/06/2024

#### 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