

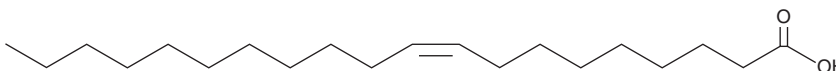
# PRODUCT INFORMATION



## Gadoleic Acid

Item No. 33186

CAS Registry No.: 29204-02-2  
Formal Name: 9Z-eicosenoic acid  
Synonyms: C20:1(9Z), cis-11-Eicosenoic Acid,  
FA 20:1  
MF:  $C_{20}H_{38}O_2$   
FW: 310.5  
Purity:  $\geq 98\%$   
Supplied as: A solution in ethanol  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 2$  years



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

### Description

Gadoleic acid is a monounsaturated fatty acid.<sup>1,2</sup> It inhibits the dsDNA binding activity of p53 when used at a concentration of 1.2 nM.<sup>1</sup> Hepatic levels of gadoleic acid are reduced in rats fed a high-fat or a high-fat high-cholesterol diet and increased in rats fed a high-cholesterol diet.<sup>2</sup>

### References

1. Iijima, H., Kasai, N., Chiku, H., *et al.* The inhibitory action of long-chain fatty acids on the DNA binding activity of p53. *Lipids* **41**(6), 521-527 (2006).
2. Serviddio, G., Bellanti, F., Villani, R., *et al.* Effects of dietary fatty acids and cholesterol excess on liver injury: A lipidomic approach. *Redox Biol.* **9**, 296-305 (2016).

#### 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, 02/20/2024

#### 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