

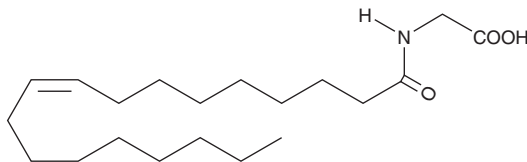
# PRODUCT INFORMATION



## N-Oleoyl Glycine

Item No. 90269

CAS Registry No.: 2601-90-3  
Formal Name: N-[(9Z)-1-oxo-9-octadecen-1-yl]-glycine  
MF:  $C_{20}H_{37}NO_3$   
FW: 339.5  
Purity:  $\geq 98\%$   
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

N-Oleoyl glycine is supplied as a crystalline solid. A stock solution may be made by dissolving the N-oleoyl glycine in the solvent of choice, which should be purged with an inert gas. N-Oleoyl glycine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N-oleoyl glycine in ethanol and DMSO is approximately 12 mg/ml and approximately 10 mg/ml in DMF.

N-Oleoyl glycine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, N-oleoyl glycine should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. N-Oleoyl glycine has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

N-Oleoyl glycine is a putative substrate for peptidyl glycine  $\alpha$ -amidating enzyme in the biosynthesis of oleoyl amide.<sup>1,2</sup>

### References

1. Merkle, D.J., Merkle, K.A., Stern, W., *et al.* Fatty acid amide biosynthesis: A possible role for peptidylglycine  $\alpha$ -amidating enzyme and acyl-coenzyme A:glycine N-acyltransferase. *Arch. Biochem. Biophys.* **330(2)**, 430-434 (1996).
2. Cravatt, B.F., Prospero-Garcia, O., Siuzdak, G., *et al.* Chemical characterization of a family of brain lipids that induce sleep. *Science* **268(5216)**, 1506-1509 (1995).

#### 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, 10/21/2022

#### 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