

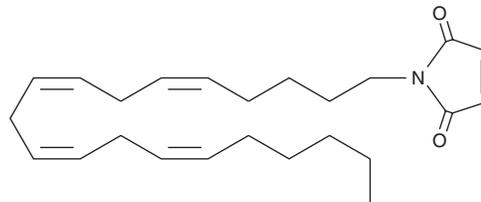
# PRODUCT INFORMATION



## N-Arachidonyl Maleimide

Item No. 10007517

**CAS Registry No.:** 876305-42-9  
**Formal Name:** eicosa-5Z,8Z,11Z,14Z-tetraenyl-1-pyrrole-2,5-dione  
**Synonym:** NAM  
**MF:** C<sub>24</sub>H<sub>35</sub>NO<sub>2</sub>  
**FW:** 369.5  
**Purity:** ≥98%  
**Supplied as:** A solution in methyl acetate  
**Storage:** -80°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

N-Arachidonyl maleimide (NAM) is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate 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 NAM in ethanol is approximately 50 mg/ml and approximately 30 mg/ml in DMSO and DMF.

### Description

2-Arachidonoyl glycerol (2-AG) is an endogenous agonist of the central cannabinoid (CB<sub>1</sub>) and peripheral cannabinoid (CB<sub>2</sub>) receptors.<sup>1-3</sup> 2-AG is present at relatively high levels in the central nervous system and is the most abundant molecular species of monoacylglycerol found in rat brain.<sup>1,4</sup> Monoacylglycerol lipase (MGL) hydrolyzes 2-AG to arachidonic acid and glycerol, thereby terminating its biological actions.<sup>5</sup> NAM is a potent, irreversible inhibitor of MGL or MGL-like activity in rat cerebellar membranes, exhibiting an IC<sub>50</sub> value of 140 nM.<sup>6</sup> Inhibition of MGL by the sulfhydryl-reactive maleimide group of NAM suggests a critical cysteine residue is present in the substrate-binding site of the enzyme.

### References

1. Stella, N., Schweitzer, P., and Piomelli, D. *Nature* **388**, 773-778 (1997).
2. Sugiura, T., Kodaka, T., Nakane, S., et al. *J. Biol. Chem.* **274**, 2794-2801 (1999).
3. Ben-Shabat, S., Frider, E., Sheskin, T., et al. *Eur. J. Pharmacol.* **353**, 23-31 (1998).
4. Kondo, S., Kondo, H., Nakane, S., et al. *FEBS Lett.* **429**, 152-156 (1998).
5. Dinh, T.P., Carpenter, D., Leslie, F.M., et al. *Proc. Natl. Acad. Sci. USA* **99**(16), 10819-10824 (2002).
6. Saario, S.M., Salo, O.M.H., Nevalainen, T., et al. *Chemistry & Biology* **12**, 649-656 (2005).

#### 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, 10/10/2023

#### 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