

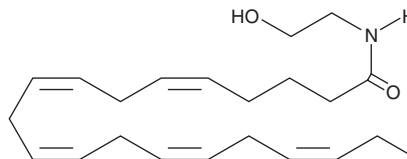
# PRODUCT INFORMATION



## Eicosapentaenoyl Ethanolamide

Item No. 10964

**CAS Registry No.:** 109001-03-8  
**Formal Name:** N-(2-hydroxyethyl)-5Z,8Z,11Z,14Z,17Z-eicosapentaenamide  
**Synonym:** EPEA  
**MF:** C<sub>22</sub>H<sub>35</sub>NO<sub>2</sub>  
**FW:** 345.5  
**Purity:** ≥98%  
**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

Eicosapentaenoyl ethanolamide (EPEA) 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 DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of EPEA in these solvents is approximately 30 mg/ml.

EPEA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of EPEA should be diluted with the aqueous buffer of choice. EPEA has a solubility of 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

EPEA is an N-acylethanolamide that inhibits dietary-restriction-induced lifespan extension in wild type and TOR pathway mutant nematodes.<sup>1</sup> Produced endogenously from eicosapentaenoic acid, EPEA serves as a metabolic signal that couples nutrient availability with growth and lifespan. EPEA also has anti-inflammatory properties, suppressing the expression of IL-6 and MCP-1 in 3T3-L1 adipocytes in response to lipopolysaccharide.<sup>2</sup>

### References

1. Lucanic, M., Held, J.M., Vantipalli, M.C., *et al.* N-acylethanolamine signalling mediates the effect of diet on lifespan in *Caenorhabditis elegans*. *Nature* **473**(7346), 226-9 (2011).
2. Balvers, M.G., Verhoeckx, K.C., Plastina, P., *et al.* Docosahexaenoic acid and eicosapentaenoic acid are converted by 3T3-L1 adipocytes to N-acyl ethanolamines with anti-inflammatory properties. *Biochim. Biophys. Acta* **1801**(10), 1107-1114 (2010).

**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, 02/27/2024

### 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