



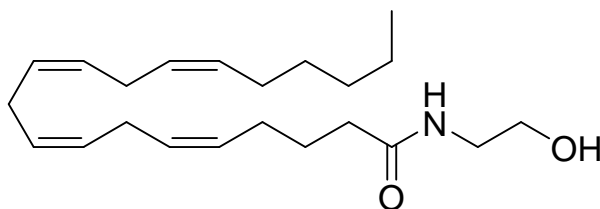
Echelon Biosciences Inc.  
675 Arapeen Drive, Suite 302  
Salt Lake City, UT 84108  
Telephone 866-588-0455  
Fax 801-588-0497  
echelon@echelon-inc.com  
www.echelon-inc.com

## Technical Data Sheet

**For research use only**  
Not intended or approved for  
diagnostic or therapeutic use.

**Product Name:** Arachidonylethanolamide (Anandamide, AEA)

**Product Number:** L-2040



**Molecular Formula:** C<sub>22</sub>H<sub>37</sub>NO<sub>2</sub>    **MW:** 347.53    **CAS:** 94421-68-8

**Solubility:** EtOH, CHCl<sub>3</sub>, Ethyl acetate: >5 mg/mL

**Storage:** -20 °C, protect from light.

**Field of Interest:** Arachidonylethanolamide is an endogenous ligand for the CB<sub>1</sub> and CB<sub>2</sub> cannabinoid receptors and vanilloid receptor. It is produced by hydrolysis of N-acyl phosphatidylethanolamine (NAPE) by NAPE-specific phospholipase D (NAPE-PLD). Fatty acid amide hydrolase (FAAH) degrades AEA to arachidonic acid and ethanolamine.

**References:** 1) W.A. Devane *et al.* "Isolation and structure of a brain constituent that binds to the cannabinoid receptor" *Science*, **1992**, 258, 1946-9. 2) J. Wang, N. Ueda "Biology of the eicocannabinoid synthesis system" *Prostaglandins and other Lipid Mediators*, **2009**, 89, 112-119.

**Hazardous Properties and Cautions:** The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. Arachidonylethanolamide is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

**Warranty and Disclaimer:** Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.