

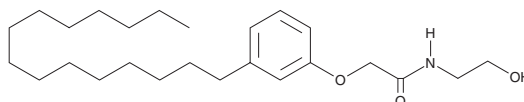
# PRODUCT INFORMATION



## PDPEA

Item No. 30956

CAS Registry No.: 861891-72-7  
Formal Name: N-(2-hydroxyethyl)-2-(3-pentadecylphenoxy)-acetamide  
MF:  $C_{25}H_{43}NO_3$   
FW: 405.6  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

PDPEA is supplied as a solid. A stock solution may be made by dissolving the PDPEA in the solvent of choice, which should be purged with an inert gas. PDPEA is soluble in organic solvents such as ethanol and DMSO. The solubility of PDPEA in these solvents is approximately 50 mM.

### Description

PDPEA is an enhancer of fatty acid amide hydrolase (FAAH) activity.<sup>1</sup> It blocks inhibition of recombinant *A. thaliana* or rat FAAH activity induced by N-acylethanolamine 12:0 (NAE 12:0) when used at a concentration of 100  $\mu\text{M}$ . PDPEA (20  $\mu\text{M}$ ) inhibits NAE 12:0-induced decreases in cotyledon size in *A. thaliana* seedlings.

### Reference

1. Faure, L., Nagarajan, S., Hwang, H., et al. Synthesis of phenoxyacyl-ethanolamides and their effects on fatty acid amide hydrolase activity. *J. Biol. Chem.* **289**(13), 9340-9351 (2014).

#### 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, 11/04/2022

#### 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