

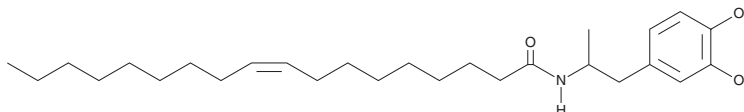
# PRODUCT INFORMATION



## N-(1-(3,4-Dihydroxyphenyl)propan-2-yl)oleamide

Item No. 9002954

**CAS Registry No.:** 1258011-97-0  
**Formal Name:** N-[2-(3,4-dihydroxyphenyl)-1-methylethyl]-9Z-octadecenamide  
**MF:** C<sub>27</sub>H<sub>45</sub>NO<sub>3</sub>  
**FW:** 431.7  
**Purity:** ≥97%  
**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

N-(1-(3,4-Dihydroxyphenyl)propan-2-yl)oleamide 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 N-(1-(3,4-dihydroxyphenyl)propan-2-yl)oleamide in these solvents is approximately 25 and 30 mg/ml, respectively.

N-(1-(3,4-Dihydroxyphenyl)propan-2-yl)oleamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of N-(1-(3,4-dihydroxyphenyl)propan-2-yl)oleamide should be diluted with the aqueous buffer of choice. N-(1-(3,4-Dihydroxyphenyl)propan-2-yl)oleamide has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method.

### Description

N-(1-(3,4-Dihydroxyphenyl)propan-2-yl)oleamide binds to the cannabinoid 1 (CB<sub>1</sub>) receptor with a K<sub>i</sub> value of 365 nM in a radioligand binding assay using rat brain homogenate.<sup>1</sup> It has an EC<sub>50</sub> value of 698 nM for the peroxisome proliferator-activated receptor α (PPARα) in a luciferase reporter assay and, in rats, it decreases food intake. It does not inhibit fatty acid amide hydrolase (FAAH).

### Reference

1. Almeida, B., Joglar, J., Rojas, M.J.L., *et al.* Synthesis of fatty acid amides of catechol metabolites that exhibit antiobesity properties. *ChemMedChem* **5**(10), 1781-1787 (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, 02/22/2024

#### 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