

PRODUCT INFORMATION



R-2 Methanandamide

Item No. 90074

CAS Registry No.: 57182-47-3

Formal Name: N-(2R-hydroxypropyl)-5Z,8Z,11Z,14Z-eicosatetraenamide

Synonym: (R)-(-)-Arachidonyl-2'-Hydroxy-1'-Propylamide

MF: $C_{23}H_{39}NO_2$

FW: 361.6

Purity: $\geq 98\%$

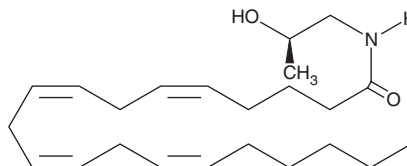
Supplied as: A 10 mg/ml solution in ethanol

Storage: -20°C

Stability: ≥ 4 years

Special Conditions: Oxygen and light sensitive

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



Laboratory Procedures

R-2 Methanandamide 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, dimethyl formamide, or acetonitrile purged with an inert gas or nitrogen can be used. Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations.

Description

R-2 Methanandamide is a cannabinoid analog with a methyl group in the (R) configuration at C-2 of the ethanolamine group. In contrast to the other methyl-anandamide analogs, R-2 methanandamide is not an amidohydrolase inhibitor and is nearly as susceptible to amide hydrolysis as AEA itself.¹

Reference

1. Abadji, V., Lin, S., Taha, G., et al. (R)-Methanandamide: A chiral novel anandamide possessing higher potency and metabolic stability. *J. Med. Chem.* **37**, 1889-1893 (1994).

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, 06/18/2026

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