

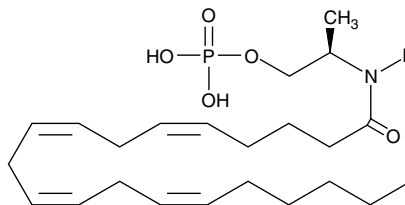
# Product Information



## R-1 Methanandamide Phosphate

Item No. 10004281

**CAS Registry No.:** 649569-33-5  
**Formal Name:** N-(2-phosphate-1R-methylethyl)-5Z,8Z,11Z,14Z-eicosatetraenamide  
**Synonyms:** R-1MAP, (R)-(+)-Arachidonyl-1'-Hydroxy-2'-Propylamide Phosphate  
**MF:** C<sub>23</sub>H<sub>40</sub>NO<sub>5</sub>P  
**FW:** 441.5  
**Purity:** ≥98%  
**Stability:** ≥1 year at -20°C  
**Supplied as:** A solution in ethanol



### Laboratory Procedures

For long term storage, we suggest that R-1 methanandamide phosphate (R-1MAP) be stored as supplied at -20°C. It should be stable for at least one year.

R-1MAP 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 R-1MAP in these solvents is at least 15 mg/ml.

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. If an organic solvent-free solution of R-1MAP is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of R-1MAP in PBS (pH 7.2) is at least 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Arachidonyl ethanolamide (AEA) was the first endogenous cannabinoid (CB) to be isolated and characterized as an agonist acting on the same receptors (CB<sub>1</sub> and CB<sub>2</sub>) as Δ<sup>9</sup>-THC.<sup>1,2</sup> Since that time, a number of related endocannabinoids have been isolated, most notably 2-arachidonoyl glycerol (2-AG).<sup>2</sup> The phosphate ester of R-1 methanandamide, R-1MAP, has been tested as a water soluble prodrug analog of AEA.<sup>3</sup> The activity of R-1MAP was essentially equivalent to that of AEA in the growth inhibition of C6 glioma cells. However, when tested for inhibition of AEA binding to isolated rat brain CB<sub>1</sub> receptors, arachidonoyl ethanolamide phosphate (AEA-P) is about 5-fold less potent as an agonist with a K<sub>i</sub> of about 200 nM.<sup>4</sup> The phosphate esters of AEA and its analogs are also structural variants of lysophosphatidic acid (LPA). However, the effects of R-1MAP on the various LPA receptors have not been tested.

### References

1. Devane, W.A., Hanus, L., Breuer, A., *et al.* Isolation and structure of a brain constituent that binds to the cannabinoid receptor. *Science* **258**, 1946-1949 (1992).
2. Felder, C.C., Briley, E.M., Axelrod, J., *et al.* Anandamide, an endogenous cannabimimetic eicosanoid, binds to the cloned human cannabinoid receptor and stimulates receptor-mediated signal transduction. *Proc. Natl. Acad. Sci. USA* **90**, 7656-7660 (1993).
3. Fowler, C.J., Jonsson, K.-O., Andersson, A., *et al.* Inhibition of C6 glioma cell proliferation by anandamide, 1-arachidonoylglycerol, and by a water soluble phosphate ester of anandamide: Variability in response and involvement of arachidonic acid. *Biochem. Pharmacol.* **66**, 757-767 (2003).
4. Sheskin, T., Hanus, L., Slager, J., *et al.* Structural requirements for binding of anandamide-type compounds to the brain cannabinoid receptor. *J. Med. Chem.* **40**, 659-667 (1997).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10004281](http://www.caymanchem.com/catalog/10004281)

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 08/27/2013

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

custserv@caymanchem.com

#### Web

[www.caymanchem.com](http://www.caymanchem.com)