

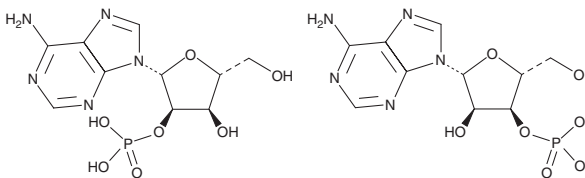
# PRODUCT INFORMATION



## Adenosine 2'(3')-monophosphate

Item No. 16062

**Formal Name:** 2'-adenylic acid, 3'-adenylic acid  
**Synonym:** 2'(3')-AMP  
**MF:** C<sub>10</sub>H<sub>14</sub>N<sub>5</sub>O<sub>7</sub>P  
**FW:** 347.2  
**Purity:** ≥95% (mixture of isomers)  
**UV/Vis.:** λ<sub>max</sub>: 259 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Adenosine 2'(3')-monophosphate is supplied as a crystalline solid. Aqueous solutions of adenosine 2'(3')-monophosphate can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of adenosine 2'(3')-monophosphate in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Adenosine 2'(3')-monophosphate is a mixture of adenosine monophosphate (AMP) derivatives with the phosphate bound at the 2'- or 3'-position as opposed to the 5'-position in AMP. Non-canonical AMP derivatives have been used to study AMP-dependent cellular processes.<sup>1,2</sup>

### References

1. Ohkubo, S., Nakanishi, H., Kimura, J., *et al.* Effects of AMP derivatives on cyclic AMP levels in NG108-15 cells. *Br. J. Pharmacol.* **129(6)**, 1244-1250 (2000).
2. Ding, J., Koellner, G., Grunert, H.P., *et al.* Crystal structure of ribonuclease T<sub>1</sub> complexed with adenosine 2'-monophosphate at 1.8-Å resolution. *J. Biol. Chem.* **266(23)**, 15128-15134 (1991).

#### 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, 05/03/2023

#### 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