

# PRODUCT INFORMATION



## N<sup>6</sup>-[(6-Aminohexyl)carbamoylmethyl]-ADP

Item No. 38402

**CAS Registry No.:** 100826-95-7  
**Formal Name:** N-[2-[(6-aminohexyl)amino]-2-oxoethyl]-adenosine 5'-(trihydrogen diphosphate)  
**Synonyms:** ACM-Adenosine 5'-diphosphate, ACM-ADP, N<sup>6</sup>-[(6-Aminohexyl)carbamoylmethyl]-Adenosine 5'-diphosphate

**MF:** C<sub>18</sub>H<sub>31</sub>N<sub>7</sub>O<sub>11</sub>P<sub>2</sub>

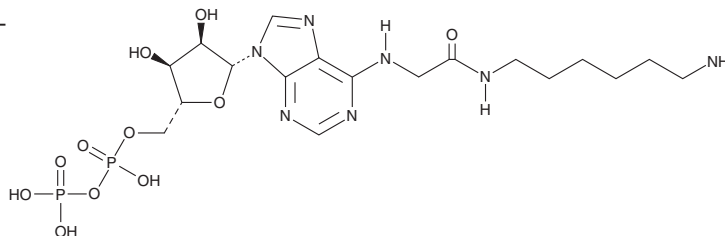
**FW:** 583.4

**Purity:** ≥70%

**Supplied as:** A crystalline solid

**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<sup>6</sup>-[(6-Aminohexyl)carbamoylmethyl]-ADP is supplied as a solid. Aqueous solutions of N<sup>6</sup>-[(6-aminohexyl)carbamoylmethyl]-ADP can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of N<sup>6</sup>-[(6-aminohexyl)carbamoylmethyl]-ADP in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

N<sup>6</sup>-[(6-Aminohexyl)carbamoylmethyl]-ADP is an alkylated derivative of the nucleotide and ATP precursor ADP (Item Nos. 16778 | 21121).<sup>1</sup> N<sup>6</sup>-[(6-Aminohexyl)carbamoyl-methyl]-ADP, bound to agarose beads, can be phosphorylated *via* acetate kinase to ATP (Item No. 14498). It has been biotinylated and conjugated to an azidonitrobenzoyl group for use as a photoaffinity probe to visualize the ATPase site of rabbit myosin.<sup>2</sup>

### References

1. Lindberg, M. and Mosbach, K. Preparation of analogues of ATP, ADP and AMP suitable for binding to matrices and the enzymic interconversion of ATP and ADP in solid phase. *Eur. J. Biochem.* **53**(2), 481-486 (1975).
2. Sutoh, K., Yamamoto, K., and Wakabayashi, T. Electron microscopic visualization of the ATPase site of myosin by photoaffinity labeling with a biotinylated photoreactive ADP analog. *Proc. Natl. Acad. Sci. USA* **83**(2), 212-216 (1986).

#### 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, 08/01/2023

#### 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