

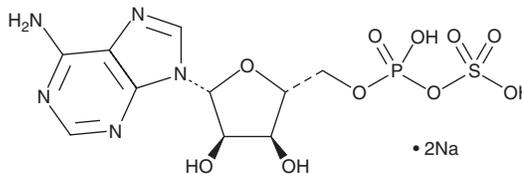
PRODUCT INFORMATION



Adenosine 5'-phosphosulfate (sodium salt)

Item No. 21226

CAS Registry No.: 102029-95-8
Formal Name: 5'-adenylic acid, sulfuric anhydride, disodium salt
Synonym: APS
MF: C₁₀H₁₄N₅O₁₀PS • 2Na
FW: 473.3
Purity: ≥95%
UV/Vis.: λ_{max}: 260 nm
Supplied as: A 10 mM aqueous solution
Storage: -20°C
Stability: ≥2 years



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

Description

Adenosine 5'-phosphosulfate is an ATP- and sulfate-competitive inhibitor of ATP sulfurylase in humans, *S. cerevisiae*, and *P. chrysogenum* ($K_i = 18, 2,500, \text{ and } 1,500 \text{ nM}$, respectively).^{1,2} Adenosine 5'-phosphosulfate also inhibits human adenosine 5'-phosphosulfate kinase ($K_i = 47.5 \mu\text{M}$) to prevent sulfation.^{2,3}

References

1. Farley, J.R., Cryns, D.F., Yang, Y.H.J., et al. Adenosine triphosphate sulfurylase from *Penicillium chrysogenum*. Steady state kinetics of the forward and reverse reactions. *J. Biol. Chem.* **251(14)**, 4389-4397 (1976).
2. Landson, E.B., Fisher, A.J., and Segel, I.H. Human 3'-phosphoadenosine 5'-phosphosulfate synthetase (isoform 1, brain): Kinetic properties of the adenosine triphosphate sulfurylase and adenosine 5'-phosphosulfate kinase domains. *Biochemistry* **43(14)**, 4656-4365 (2004).
3. Mueller, J.W. and Shafqat, N. Adenosine-5'-phosphosulfate--a multifaceted modulator of bifunctional 3'-phospho-adenosine-5'-phosphosulfate synthases and related enzymes. *FEBS J.* **280(13)**, 3050-3057 (2013).

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, 04/25/2018

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