

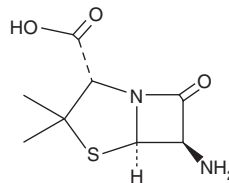
PRODUCT INFORMATION



6-Aminopenicillanic Acid

Item No. 34494

CAS Registry No.: 551-16-6
Formal Name: (2S,5R,6R)-6-amino-3,3-dimethyl-7-oxo-4-thia-1-azabicyclo[3.2.0]heptane-2-carboxylic acid
Synonyms: (+)-6-Aminopenicillanic Acid, 6 β -Aminopenicillanic Acid, 6-APA, NSC 50071
MF: C₈H₁₂N₂O₃S
FW: 216.3
Purity: $\geq 90\%$
Supplied as: A 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

6-Aminopenicillanic acid (6-APA) is supplied as a solid. Aqueous solutions of 6-APA can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 6-APA in PBS (pH 7.2) is approximately 1 mg/ml (sonicated). We do not recommend storing the aqueous solution for more than one day.

Description

6-APA is an intermediate in the semisynthetic synthesis of penicillin antibiotics.¹

Reference

1. Su, M., Sun, H., Zhao, Y., *et al.* Degradation kinetics and mechanism of a β -lactam antibiotic intermediate, 6-aminopenicillanic acid, in a new integrated production process. *J. Pharm. Sci.* **105**(1), 139-146 (2016).

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, 11/22/2022

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