

PRODUCT INFORMATION



SAICAR (hydrate)

Item No. 14087

Formal Name: N-[[5-amino-1-(5-O-phosphono-β-D-ribofuranosyl)-1H-imidazol-4-yl]carbonyl]-L-aspartic acid, hydrate

Synonyms: Succino-AICAR, Succinyl-5-aminoimidazole-4-carboxamide-1-ribose-5'-phosphate

MF: C₁₃H₁₉N₄O₁₂P • XH₂O

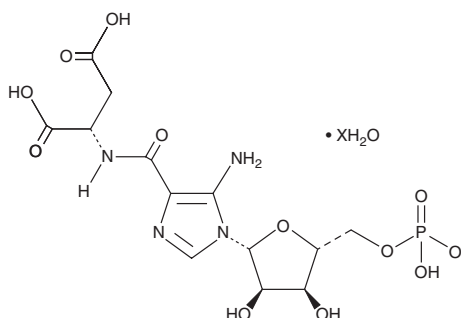
FW: 454.3

Purity: ≥98%

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

SAICAR (hydrate) is supplied as a solid. A stock solution may be made by dissolving the SAICAR (hydrate) in the solvent of choice, which should be purged with an inert gas. SAICAR (hydrate) is slightly soluble in methanol.

SAICAR (hydrate) is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

SAICAR is an intermediate in the *de novo* purine nucleotide synthesis pathway. It is reported to stimulate pyruvate kinase isoform M2 and to promote cancer cell survival in conditions of limited glucose availability.¹

Reference

1. Keller, K.E., Tan, I.S., and Lee, Y.-S. SAICAR stimulates pyruvate kinase isoform M2 and promotes cancer cell survival in glucose-limited conditions. *Science* **338**(6110), 1069-1072 (2012).

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, 12/01/2022

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