

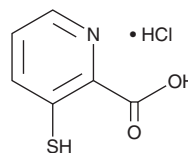
PRODUCT INFORMATION



3-Mercaptopicolinic Acid (hydrochloride)

Item No. 20895

CAS Registry No.: 320386-54-7
Formal Name: 3-mercapto-2-pyridinecarboxylic acid, monohydrochloride
Synonym: SKF 34288
MF: C₆H₅NO₂S • HCl
FW: 191.6
Purity: ≥95%
UV/Vis.: λ_{max}: 209, 251, 316, 431 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

3-Mercaptopicolinic acid (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of 3-mercapto-2-pyridinecarboxylic acid (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 3-mercapto-2-pyridinecarboxylic acid (hydrochloride) in PBS (pH 7.2) is approximately 0.2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

3-Mercaptopicolinic acid is an inhibitor of gluconeogenesis. In perfused rat livers, gluconeogenesis was inhibited when lactate, pyruvate, or alanine served as substrate, but not with fructose, suggesting phosphoenolpyruvate carboxykinase as the site of inhibition.¹⁻³

References

1. DiTullio, N.W., Berkoff, C.E., Blank, B., *et al.* 3-Mercaptopicolinic acid, an inhibitor of gluconeogenesis. *Biochem. J.* **138(3)**, 387-394 (1974).
2. Robinson, B.H. and Oei, J. 3-Mercaptopicolinic acid, a preferential inhibitor of the cytosolic phosphoenolpyruvate carboxykinase. *FEBS Lett.* **58(1)**, 12-15 (1975).
3. Hidalgo, J., Latorre, P., Carrodeguas, J.A., *et al.* Inhibition of pig phosphoenolpyruvate carboxykinase isoenzymes by 3-mercapto-2-pyridinecarboxylic acid and novel inhibitors. *PLoS One* **11(7)**, e0159002 (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, 11/04/2022

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