

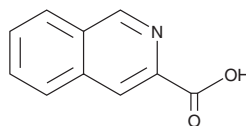
PRODUCT INFORMATION



Isoquinoline-3-carboxylic Acid

Item No. 19194

CAS Registry No.: 6624-49-3
Formal Name: 3-isoquinolinecarboxylic acid
Synonyms: 3-Carboxyisoquinoline,
3-Isoquinaldic Acid, NSC 53385
MF: C₁₀H₇NO₂
FW: 173.2
Purity: ≥98%
UV/Vis.: λ_{max}: 229 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

Isoquinoline-3-carboxylic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the isoquinoline-3-carboxylic acid in the solvent of choice, which should be purged with an inert gas. Isoquinoline-3-carboxylic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of isoquinoline-3-carboxylic acid in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Isoquinoline-3-carboxylic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, isoquinoline-3-carboxylic acid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Isoquinoline-3-carboxylic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Isoquinoline-3-carboxylic acid is a synthetic intermediate that has been used to design novel antitumor compounds.¹

Reference

1. Gao, F., Liu, H., Li, L., *et al.* Design, synthesis, and testing of an isoquinoline-3-carboxylic-based novel anti-tumor lead. *Bioorg. Med. Chem. Lett.* **25(20)**, 4434-4436 (2015).

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, 11/10/2022

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