

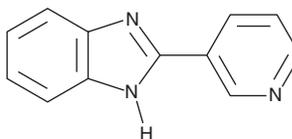
PRODUCT INFORMATION



2-(3-pyridyl)-Benzimidazole

Item No. 13054

CAS Registry No.: 1137-67-3
Formal Name: 2-(3-pyridinyl)-1H-benzimidazole
Synonyms: G 514, NSC 62610
MF: C₁₂H₉N₃
FW: 195.2
Purity: ≥98%
UV/Vis.: λ_{max}: 307 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

2-(3-pyridyl)-Benzimidazole is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-(3-pyridyl)-benzimidazole in an organic solvent purged with an inert gas. 2-(3-pyridyl)-Benzimidazole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2-(3-pyridyl)-benzimidazole in ethanol is approximately 15 mg/ml and approximately 10 mg/ml in DMSO and DMF.

2-(3-pyridyl)-Benzimidazole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2-(3-pyridyl)-benzimidazole should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 2-(3-pyridyl)-Benzimidazole has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

2-(3-pyridyl)-Benzimidazole is a synthetic intermediate useful for pharmaceutical synthesis.

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, 12/13/2022

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