

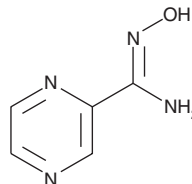
PRODUCT INFORMATION



Pyrazine-2-amidoxime

Item No. 11059

CAS Registry No.: 51285-05-3
Formal Name: N-hydroxy-2-pyrazinecarboximidamide
MF: C₅H₆N₄O
FW: 138.1
Purity: ≥97%
UV/Vis.: λ_{max}: 206, 251, 302 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

Pyrazine-2-amidoxime is supplied as a crystalline solid. A stock solution may be made by dissolving the pyrazine-2-amidoxime in an organic solvent purged with an inert gas. Pyrazine-2-amidoxime is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of pyrazine-2-amidoxime in ethanol is approximately 0.2 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of pyrazine-2-amidoxime can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of pyrazine-2-amidoxime in PBS, pH 7.2, is approximately 0.1 mg/ml.

Description

Pyrazine-2-amidoxime 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, 12/12/2022

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