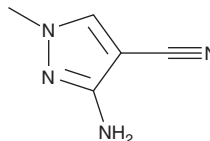


PRODUCT INFORMATION

1-Methyl-3-amino-4-cyanopyrazole

Item No. 42224

CAS Registry No.: 21230-50-2
 Formal Name: 3-amino-1-methyl-1H-pyrazole-4-carbonitrile
 MF: $C_5H_6N_4$
 FW: 122.1
 Purity: $\geq 98\%$
 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

1-Methyl-3-amino-4-cyanopyrazole is supplied as a solid. A stock solution may be made by dissolving the 1-methyl-3-amino-4-cyanopyrazole in the solvent of choice, which should be purged with an inert gas. 1-Methyl-3-amino-4-cyanopyrazole is soluble (≥ 10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

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 1-methyl-3-amino-4-cyanopyrazole can be prepared by directly dissolving the solid in aqueous buffers. 1-Methyl-3-amino-4-cyanopyrazole is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

1-Methyl-3-amino-4-cyanopyrazole is a precursor in the synthesis of adenosine A_3 receptor antagonists with anticancer activity *in vitro*.¹

Reference

1. Taliani, S., La Motta, C., Mugnaini, L., *et al.* Novel N^2 -substituted pyrazolo[3,4-d]pyrimidine adenosine A_3 receptor antagonists: Inhibition of A_3 -mediated human glioblastoma cell proliferation. *J. Med. Chem.* **53**(10), 3954-3963 (2010).

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, 11/07/2024

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