

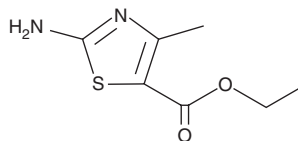
PRODUCT INFORMATION



ethyl-2-amino-4-methyl-Thiazole-5-Carboxylate

Item No. 13059

CAS Registry No.: 7210-76-6
Formal Name: 2-amino-4-methyl-5-thiazolecarboxylic acid, ethyl ester
Synonyms: EMTC, NSC 400231
MF: C₇H₁₀N₂O₂S
FW: 186.2
Purity: ≥98%
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

ethyl-2-amino-4-methyl-Thiazole-5-carboxylate is supplied as a crystalline solid. A stock solution may be made by dissolving the ethyl-2-amino-4-methyl-thiazole-5-carboxylate in the solvent of choice, which should be purged with an inert gas. ethyl-2-amino-4-methyl-Thiazole-5-carboxylate is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

ethyl-2-amino-4-methyl-Thiazole-5-carboxylate 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