

# PRODUCT INFORMATION



## 2-amino-5-methyl-4-phenyl-Thiazole

Catalog No. 13061

CAS Registry No.: 30709-67-2

Formal Name: 5-methyl-4-phenyl-2-thiazolamine

Synonym: NSC 54435

MF:  $C_{10}H_{10}N_2S$

FW: 190.3

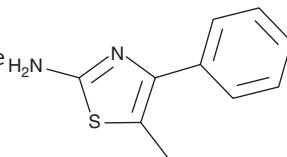
Purity:  $\geq 98\%$

UV/Vis.:  $\lambda_{max}$ : 224, 260 nm

Supplied as: A crystalline solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

2-amino-5-methyl-4-phenyl-Thiazole is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-amino-5-methyl-4-phenyl-thiazole in the organic solvent of choice, which should be purged with an inert gas. 2-amino-5-methyl-4-phenyl-Thiazole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2-amino-5-methyl-4-phenyl-thiazole in ethanol is approximately 12.5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

2-amino-5-methyl-4-phenyl-Thiazole is sparingly soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. For biological experiments, we suggest that organic solvent-free aqueous solutions of 2-amino-5-methyl-4-phenyl-thiazole be prepared by directly dissolving the 2-amino-5-methyl-4-phenyl-thiazole compound in aqueous buffers. The solubility of 2-amino-5-methyl-4-phenyl-thiazole in PBS, pH 7.2, is approximately 0.15 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

2-amino-5-methyl-4-phenyl-Thiazole 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