

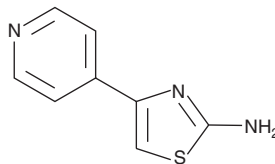
# PRODUCT INFORMATION



## 2-amino-4-(4-pyridyl)-Thiazole

Item No. 11058

CAS Registry No.: 30235-28-0  
Formal Name: 4-(4-pyridinyl)-2-thiazolamine  
MF:  $C_8H_7N_3S$   
FW: 177.2  
Purity:  $\geq 97\%$   
UV/Vis.:  $\lambda_{max}$ : 233, 314 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-4-(4-pyridyl)-Thiazole is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-amino-4-(4-pyridyl)-thiazole in the organic solvent of choice, which should be purged with an inert gas. 2-amino-4-(4-pyridyl)-Thiazole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2-amino-4-(4-pyridyl)-Thiazole in these solvents is approximately 0.14, 10, and 5 mg/ml, respectively.

2-amino-4-(4-pyridyl)-Thiazole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2-amino-4-(4-pyridyl)-thiazole should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 2-amino-4-(4-pyridyl)-Thiazole has a solubility of approximately 0.1 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

2-amino-4-(4-pyridyl)-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, 12/12/2022

#### 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