

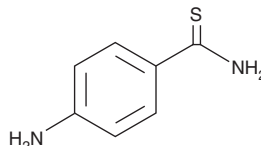
PRODUCT INFORMATION



4-amino Benzthioamide

Item No. 11063

CAS Registry No.: 4714-67-4
Synonym: 4-Aminobenzenecarbothioamide
MF: $C_7H_8N_2S$
FW: 152.2
Purity: $\geq 97\%$
UV/Vis.: λ_{max} : 338 nm
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

4-amino Benzthioamide is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-amino benzthioamide in the solvent of choice, which should be purged with an inert gas. 4-amino Benzthioamide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 4-amino benzthioamide in ethanol is approximately 0.15 mg/ml and approximately 10 mg/ml in DMSO and DMF.

4-amino Benzthioamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4-amino benzthioamide should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 4-amino Benzthioamide has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

4-amino Benzthioamide 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, 10/25/2022

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