

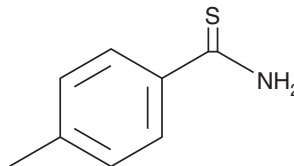
PRODUCT INFORMATION



thio-*p*-Toluamide

Item No. 13229

CAS Registry No.: 2362-62-1
Formal Name: 4-methyl-benzenecarbothioamide
Synonym: 4-Methylbenzenethioamide
MF: C₈H₉NS
FW: 151.2
Purity: ≥98%
UV/Vis.: λ_{max}: 255, 301 nm
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

thio-*p*-Toluamide is supplied as a crystalline solid. A stock solution may be made by dissolving the thio-*p*-toluamide in the solvent of choice, which should be purged with an inert gas. thio-*p*-Toluamide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of thio-*p*-toluamide in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

thio-*p*-Toluamide is a cyclic thioamide building block used in drug discovery chemistry.

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, 11/08/2022

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