

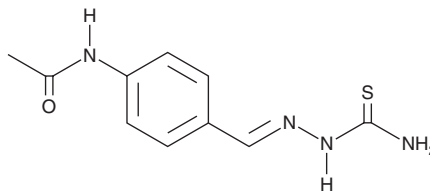
PRODUCT INFORMATION



Thioacetazone

Item No. 34589

CAS Registry No.: 104-06-3
Formal Name: N-[4-[[2-(aminothioxomethyl)hydrazinylidene]methyl]phenyl]-acetamide
Synonyms: Thiacetazone, NSC 3550, SQ 2,321
MF: C₁₀H₁₂N₄OS
FW: 236.3
Purity: ≥98%
UV/Vis.: λ_{max}: 330 nm
Supplied as: A 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

Thioacetazone is supplied as a solid. A stock solution may be made by dissolving the thioacetazone in the solvent of choice, which should be purged with an inert gas. Thioacetazone is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of thioacetazone in these solvents is approximately 30 mg/ml.

Thioacetazone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, thioacetazone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Thioacetazone has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Thioacetazone is an antitubercular agent.¹ It is active against *M. tuberculosis* H37Rv, *M. kansasii* Bostrom, and *M. avium* serotype 2 (MICs = 0.3, 0.6, and 20 µg/ml, respectively) but not *M. simiae* N29 or *M. intracellulare* serotype 16 (MICs = >40 µg/ml for both). Formulations containing thioacetazone were previously used in the treatment of tuberculosis.

Reference

1. Collins, F.M., Klayman, D.L., and Morrison, N.E. Activity of 2-acetylpyridine and 2-acetylquinoline thiosemicarbazones tested *in vitro* in combination with other antituberculous drugs. *Am. Rev. Respir. Dis.* **125**(1), 58-60 (1982).

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, 12/06/2022

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