

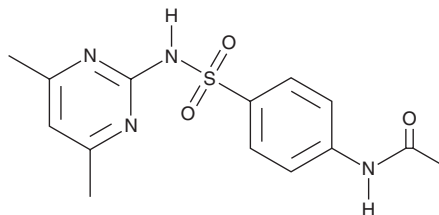
PRODUCT INFORMATION



N-acetyl Sulfamethazine

Item No. 42996

CAS Registry No.: 100-90-3
Formal Name: N-[4-[[[4,6-dimethyl-2-pyrimidinyl]amino]sulfonyl]phenyl]-acetamide
MF: $C_{14}H_{16}N_4O_3S$
FW: 320.4
Purity: $\geq 98\%$
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

N-acetyl Sulfamethazine is supplied as a solid. A stock solution may be made by dissolving the N-acetyl sulfamethazine in the solvent of choice, which should be purged with an inert gas. N-acetyl Sulfamethazine is slightly soluble (0.1-1 mg/ml) in DMSO and acetonitrile.

Description

N-acetyl Sulfamethazine is an inactive metabolite of the sulfonamide antibiotic sulfamethazine (Item No. 20976).¹

Reference

1. López-Serna, R., Petrović, M., and Barceló, D. Direct analysis of pharmaceuticals, their metabolites and transformation products in environmental waters using on-line TurboFlow™ chromatography-liquid chromatography-tandem mass spectrometry. *J. Chromatogr. A.* **1252**, 115-129 (2012).

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, 03/18/2025

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