

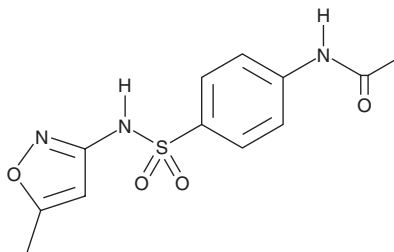
PRODUCT INFORMATION



N-acetyl Sulfamethoxazole

Item No. 17247

CAS Registry No.: 21312-10-7
Formal Name: N-[4-[[[(5-methyl-3-isoxazolyl)amino]sulfonyl]phenyl]-acetamide
Synonym: STX 606
MF: C₁₂H₁₃N₃O₄S
FW: 295.3
Purity: ≥95%
UV/Vis.: λ_{max}: 208, 265 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

N-acetyl Sulfamethoxazole is supplied as a crystalline solid. A stock solution may be made by dissolving the N-acetyl sulfamethoxazole in the solvent of choice, which should be purged with an inert gas. N-acetyl Sulfamethoxazole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N-acetyl sulfamethoxazole in ethanol is approximately 0.25 mg/ml and approximately 50 mg/ml in DMSO and DMF.

N-acetyl Sulfamethoxazole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, N-acetyl sulfamethoxazole should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. N-acetyl Sulfamethoxazole has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

N-acetyl Sulfamethoxazole is a metabolite of sulfamethoxazole, an antibiotic with antiviral activity, that can be detected in urine.^{1,2}

References

1. Bodro, M. and Paterson, D.L. Has the time come for routine trimethoprim-sulfamethoxazole prophylaxis in patients taking biologic therapies? *Clin. Infect. Dis.* **56(11)**, 1621-1628 (2013).
2. van der Ven, A.J.A.M., Vree, T.B., van Ewijk-Beneken Kolmer, E.W.J., et al. Urinary recovery and kinetics of sulphamethoxazole and its metabolites in HIV-seropositive patients and healthy volunteers after a single oral dose of sulphamethoxazole. *Br. J. Clin. Pharmacol.* **39(6)**, 621-625 (1995).

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, 12/09/2022

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