

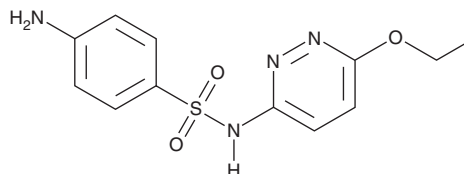
PRODUCT INFORMATION



Sulfaethoxypyridazine

Item No. 41698

CAS Registry No.: 963-14-4
Formal Name: 4-amino-N-(6-ethoxy-3-pyridazinyl)-benzenesulfonamide
MF: C₁₂H₁₄N₄O₃S
FW: 294.3
Purity: ≥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

Sulfaethoxypyridazine is supplied as a solid. A stock solution may be made by dissolving the sulfaethoxypyridazine in the solvent of choice, which should be purged with an inert gas. Sulfaethoxypyridazine is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Sulfaethoxypyridazine is a sulfonamide antibiotic.¹ It is active against several strains of *E. coli* isolated from necropsied animals or poultry when used at concentrations of 0.25 or 1 mg/ml. Prolonged administration of sulfaethoxypyridazine (150 mg/kg for 13-27 weeks) induces the formation of cataracts in dogs.² Formulations containing sulfaethoxypyridazine have been used in the treatment of respiratory infections, foot rot, calf scours, and mastitis- or metritis-associated septicemia in cattle.

References

1. Glantz, P.J. *In vitro* sensitivity of *Escherichia coli* to sulfonamides. *Cornell Vet.* **55**, 9-18 (1965).
2. Ribelin, W.E., Owen, G., Rubin, L.F., *et al.* Development of cataracts in dogs and rats from prolonged feeding of sulfaethoxypyridazine. *Toxicol. Appl. Pharmacol.* **10(3)**, 557-567 (1967).

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, 10/07/2024

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