

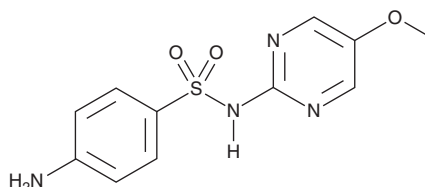
PRODUCT INFORMATION



Sulfameter

Item No. 27630

CAS Registry No.: 651-06-9
Formal Name: 4-amino-N-(5-methoxy-2-pyrimidinyl)-benzenesulfonamide
Synonyms: Bayrena, NSC 683528, Sulfamethoxydiazine
MF: C₁₁H₁₂N₄O₃S
FW: 280.3
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 230, 272 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

Sulfameter is supplied as a crystalline solid. A stock solution may be made by dissolving the sulfameter in the solvent of choice, which should be purged with an inert gas. Sulfameter is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of sulfameter in these solvents is approximately 20 and 25 mg/ml, respectively.

Sulfameter is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, sulfameter should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Sulfameter has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Sulfameter is a long-acting sulfonamide antibiotic.¹ *In vitro*, it is active against the bacteria *D. pneumoniae*, *K. pneumoniae*, *E. coli*, *N. meningitidis*, *P. vulgaris*, *S. aureus*, *S. haemolyticus*, and *S. typhimurium* in a disc assay.² Sulfameter is efficacious against lethal *K. pneumoniae*, *E. coli*, *P. vulgaris*, *D. pneumoniae*, *P. aeruginosa*, and *S. typhimurium* infection in a mouse model of septicemia with 50% curative dose (CD₅₀) values of 1.6, 5.2, 5.4, 120.5, 19.1, and 88 mg/kg, respectively. Formulations containing sulfameter have previously been used in the treatment of urinary tract infections.

References

1. Gibson, G.R. Clinical evaluation of sulphamethoxydiazine. *Med. J. Aust.* **2**, 172-174 (1964).
2. Böhni, E., Fust, B., Rieder, J., et al. Comparative toxicological, chemotherapeutic and pharmacokinetic studies with sulphormethoxine and other sulphonamides in animals and man. *Chemotherapy* **14**(4), 195-226 (1969).

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, 12/02/2022

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