PRODUCT INFORMATION



Sulfamethazine-d₄

Item No. 36575

CAS Registry No.:	1020719-82-7	
Formal Name:	4-amino-N-(4,6-dimethyl-2-pyrimidinyl)-	
	benzene-2,3,5,6-d₄-sulfonamide	Ĥ
Synonyms:	Sulfadimethyldiazine- d_4 , Sulfadimidine- d_4	
MF:	$C_{12}H_{10}D_4N_4O_2S$	
FW:	282.4	
Chemical Purity:	≥98% Sulfamethazine	
Deuterium		Han
Incorporation:	≥99% deuterated forms (d ₁ -d ₄); ≤1% d ₀	
Supplied as:	A solid	D
Storage:	-20°C	
Stability:	≥4 years	

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Sulfamethazine- d_4 is intended for use as an internal standard for the quantification of sulfamethazine (Item No. 20976) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Sulfamethazine- d_4 is supplied as a solid. A stock solution may be made by dissolving the sulfamethazine- d_4 in the solvent of choice, which should be purged with an inert gas. Sulfamethazine-d₄ is slightly soluble in DMSO and methanol (warmed).

Description

Sulfamethazine is a sulfonamide antibiotic.^{1,2} It inhibits dihydropteroate synthase (DHPS; IC_{50} = 5.7 μ M for the T. gondii enzyme). Sulfamethazine is active against A. pleuropneumoniae (MIC = 32 μ g/ml) and enhances the antibacterial activity of trimethoprim (Item No. 16473) against E. coli.^{3,4} It has been detected in environmental water samples.^{5,6} Formulations containing sulfamethazine have been used in the treatment of bacterial infections in livestock.

References

- 1. Allegra, C.J., Boarman, D., Kovacs, J.A., et al. Interaction of sulfonamide and sulfone compounds with Toxoplasma gondii dihydropteroate synthase. J. Clin. Invest. 85(2), 371-379 (1990).
- 2. Salmon, S.A., Watts, J.L., Case, C.A., et al. Comparison of MICs of ceftiofur and other antimicrobial agents against bacterial pathogens of swine from the United States, Canada, and Denmark. J. Clin. Microbiol. 33(9), 2435-2444 (1995).
- 3. Mengelers, M.J., Hougee, P.E., Janssen, L.H., et al. Structure-activity relationships between antibacterial activities and physicochemical properties of sulfonamides. J. Vet. Pharmacol. Ther. 20(4), 276-283 (1997).
- 4. Peng, F.-J., Ying, G.-G., Liu, Y.-S., et al. Joint antibacterial activity of soil-adsorbed antibiotics trimethoprim and sulfamethazine. Sci. Total Environ. 506-507, 58-65 (2015).
- 5. Washington, M.T., Moorman, T.B., Soupir, M.L., et al. Monitoring tylosin and sulfamethazine in a tiledrained agricultural watershed using polar organic chemical integrative sampler (POCIS). Sci. Total. Environ. 612, 358-367 (2017).
- 6. 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. Chomatogr. A. 1252, 115-129 (2012).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

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

uyer 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, 12/20/2022

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