

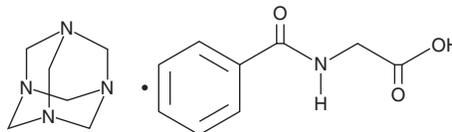
PRODUCT INFORMATION



Methenamine Hippurate

Item No. 24004

CAS Registry No.: 5714-73-8
Formal Name: N-benzoyl-glycine, compd. with 1,3,5,7-tetraazatricyclo[3.3.1.1^{3,7}]decane
MF: C₆H₁₂N₄ • C₉H₉NO₃
FW: 319.4
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

Methenamine hippurate is supplied as a solid. A stock solution may be made by dissolving the methenamine hippurate in the solvent of choice, which should be purged with an inert gas. Methenamine hippurate is soluble in organic solvents such as methanol and DMSO.

Description

Methenamine hippurate is a broad-spectrum urinary antiseptic agent.¹ Under acidic conditions, methenamine hippurate is hydrolyzed to formaldehyde and ammonia, which exerts an antibacterial effect. It is effective against a wide range of bacteria but lacks activity against *Proteus* and other bacteria that increase urine alkalinity. Formulations containing methenamine hippurate have been used to treat urinary tract infections.

Reference

1. Papich, M.G. Methenamine. *Handbook of veterinary drugs: Small and large animal*, 4th ed, Saunders, 505-506 (2016).

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, 11/16/2022

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