

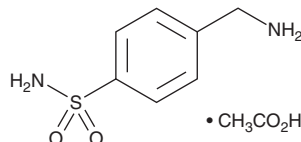
PRODUCT INFORMATION



Mafenide (acetate)

Item No. 44225

CAS Registry No.: 13009-99-9
Formal Name: 4-(aminomethyl)-benzenesulfonamide, monoacetate
Synonyms: *p*-Aminomethylbenzenesulfonamide, 4-Homosulfanilamide, Homosulfamine, Marfanil, Mesudin
MF: $C_7H_{10}N_2O_2S \cdot C_2H_4O_2$
FW: 246.3
Purity: $\geq 98\%$
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

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of mafenide (acetate) can be prepared by directly dissolving the solid in aqueous buffers. Mafenide (acetate) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Mafenide is a sulfonamide antibiotic.^{1,2} It is active against clinical isolates of *S. pyogenes*, methicillin-susceptible *S. aureus* (MSSA), methicillin-resistant *S. aureus* (MRSA), *Enterococcus*, Enterobacteriaceae, and Gram-negative bacilli from burn patients in an agar well diffusion assay (mean zone of inhibition = 24-37 mm) but not in a broth dilution assay with MIC values ranging from 250 to greater than 5,000 $\mu\text{g/ml}$.¹ Mafenide is also active against clinical isolates of *K. pneumoniae* that produce extended spectrum β -lactamase (ESBL), *P. aeruginosa*, and *A. baumannii-calcoaceticus* from burn patients in an agar well diffusion assay (mean zones of inhibition = 23.5, 28.9, and 25.8 mm, respectively) but not in a broth dilution assay (mean MICs = 1,024 $\mu\text{g/ml}$ for all).² It decreases mortality in a rat model of burn wounds seeded with rat virulent *P. aeruginosa*.³ Mafenide also inhibits human carbonic anhydrase I (CAI) and CAII (K_i s = 41.91 and 0.612 μM , respectively).⁴ Formulations containing mafenide have been used in the treatment of bacterial infections associated with burn wounds.

References

1. Rodgers, G.L., Mortensen, J.E., Fisher, M.C., et al. *J. Burn Care Rehabil.* **18**(5), 406-410 (1997).
2. Glasser, J.S., Guymon, C.H., Mende, K., et al. *Burns* **36**(8), 1172-1184 (2010).
3. Fox, C.L., Jr., Sampath, A.C., and Stanford, J.W. *Arch. Surg.* **101**(4), 508-512 (1970).
4. Fidan, İ., Salmas, R.E., Arslan, M., et al. *Bioorg. Med. Chem.* **23**(23), 7353-7358 (2015).

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, 10/14/2025

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