

PRODUCT INFORMATION

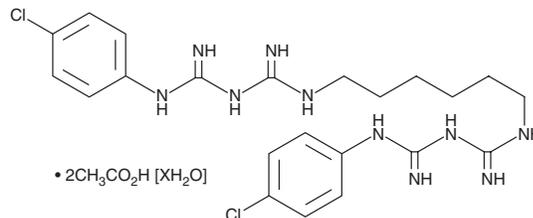


Chlorhexidine (acetate hydrate)

Item No. 26924

CAS Registry No.: 206986-79-0
Formal Name: N,N''-bis(4-chlorophenyl)-3,12-diimino-2,4,11,13-tetraazatetradecanediiimidamide, diacetate, hydrate

Synonym: CHX
MF: C₂₂H₃₀Cl₂N₁₀ • 2C₂H₄O₂ [XH₂O]
FW: 625.6
Purity: ≥98%
UV/Vis.: λ_{max}: 261 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

Chlorhexidine (acetate hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the chlorhexidine (acetate hydrate) in the solvent of choice, which should be purged with an inert gas. Chlorhexidine (acetate hydrate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of chlorhexidine (acetate hydrate) in ethanol is approximately 10 mg/ml and approximately 15 mg/ml in DMSO and DMF.

Chlorhexidine (acetate hydrate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, chlorhexidine (acetate hydrate) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Chlorhexidine (acetate hydrate) has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Chlorhexidine is a bis(biguanide) antimicrobial disinfectant and antiseptic agent.¹ It inhibits growth of clinical methicillin-resistant *S. aureus* (MRSA) isolates (MIC₉₀ = 4 µg/ml).² It is also active against canine isolates of MRSA, methicillin-susceptible *S. aureus* (MSSA), methicillin-resistant *S. pseudintermedius* (MRSP), and methicillin-susceptible *S. pseudintermedius* (MSSP; MIC₉₀s = 4, 2, 2, and 1 mg/L, respectively).³ Chlorhexidine inhibits growth of *E. faecium* strains (MICs = 1.2-19.6 µg/ml) and *C. albicans* (MIC = 5.15 µg/ml).^{4,5} It generates cations that bind to and destabilize the bacterial cell wall to induce death.⁶ Chlorhexidine also completely inhibits matrix metalloproteinase-2 (MMP-2) and MMP-9 when used at concentrations of 0.0001 and 0.002%, respectively, in a gelatin degradation assay.⁷ Formulations containing chlorhexidine have been used in antiseptic wound dressings, mouthwash, and toothpaste.

References

1. Karpiński, T.M. and Szkaradkiewicz, A.K. *Eur. Rev. Med. Pharmacol. Sci.* **19(7)**, 1321-1326 (2015).
2. McDanel, J.S., Murphy, C.R., Diekema, D.J., et al. *Antimicrob. Agents Chemother.* **57(1)**, 552-558 (2013).
3. Clark, S.M., Loeffler, A., and Bond, R. *J. Antimicrob. Chemother.* **70(7)**, 2048-2052 (2015).
4. Bhardwaj, P., Hans, A., Ruikar, K., et al. *Antimicrob. Agents Chemother.* **62(1)**, e01235-17 (2017).
5. Scheibler, E., da Silva, R.M., Leite, C.E., et al. *Arch. Oral Biol.* **89**, 70-76 (2018).
6. Barrett-Bee, K., Newbould, L., and Edwards, S. *FEMS Microbiol. Lett.* **119(1-2)**, 249-253 (1994).
7. Gendron, R., Grenier, D., Sorsa, T., et al. *Clin. Diagn. Lab. Immunol.* **6(3)**, 437-439 (1999).

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, 11/18/2022

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