

PRODUCT INFORMATION



Thimerosal

Item No. 33432

CAS Registry No.: 54-64-8

Formal Name: ethyl[2-(mercapto-κS)benzoato(2-)-κO]-mercurate(1-), monosodium

Synonyms: Merthiolate, NSC 4794

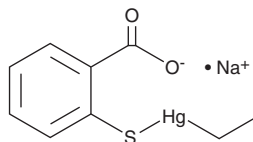
MF: $C_9H_9HgO_2S \cdot Na$

FW: 404.8

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

Thimerosal is supplied as a crystalline solid. A stock solution may be made by dissolving the thimerosal in the solvent of choice, which should be purged with an inert gas. Thimerosal is soluble in organic solvents such as ethanol. Thimerosal is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Thimerosal is a mercurial antimicrobial agent.^{1,2} It is bacteriostatic against hemolytic streptococci and fungistatic against *H. capsulatum*, *B. dermatitidis*, and *S. schenckii*. Thimerosal (1-250 μM) induces apoptosis in HCN-1A human cortical neurons and various neuroblastoma cells.^{3,4} *In vivo*, thimerosal (80 mg/kg, i.p.) induces footpad swelling and dorsal skin contact dermatitis in a mouse model of pseudo-allergic reaction.⁵ Formulations containing thimerosal have been used as antiseptics and preservatives in vaccines.

References

1. Morton, H.E., North, L.L., Jr., and Engley, F.B., Jr. The bacteriostatic and bactericidal actions of some mercurial compounds on hemolytic streptococci. *J. Am. Med. Assoc.* **136**(1), 37-41 (1948).
2. Deighton, F.J., Hall, N.K., and Larsh, H.W. Merthiolate treatment of pathogenic fungi. *J. Clin. Microbiol.* **10**(2), 144-146 (1979).
3. Baskin, D.S., Ngo, H., and Didenko, V.V. Thimerosal induces DNA breaks, caspase-3 activation, membrane damage, and cell death in cultured human neurons and fibroblasts. *Toxicol. Sci.* **74**(2), 361-368 (2003).
4. Dórea, J.G. Integrating experimental (in vitro and in vivo) neurotoxicity studies of low-dose thimerosal relevant to vaccines. *Neurochem. Res.* **36**(3), 927-938 (2011).
5. Peng, B., Che, D., Hao, Y., et al. Thimerosal induces skin pseudo-allergic reaction via Mas-related G-protein coupled receptor B2. *J. Dermatol. Sci.* **95**(3), 99-106 (2019).

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, 12/16/2022

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