

Thimerosal

Antiseptic and antifungal agent

Modulates thiol groups in biological systems, especially in those involving cAMP, Ca²⁺ and inositol phosphate. Causes a release of calcium from intracellular stores in many cells types. Has toxic effects on several cell types. Induces apoptosis *in vitro*.

Citations: 10

[View Online »](#)

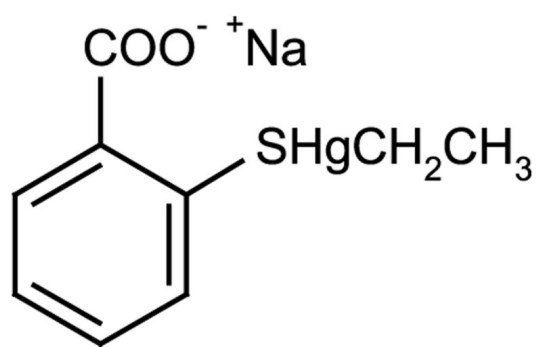
Ordering Information

[Order Online »](#)

ALX-400-013-G005	5g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ethyl mercurithiosalicylic acid . Na, Mercuriothiolate
Appearance	White to off-white powder.
CAS	54-64-8
Formula	$C_9H_9HgO_2SNa$
MI	14: 9317
MW	404.8
Purity	≥98% (HPLC)
Solubility	Soluble in water (600mg/ml) or methanol.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)