

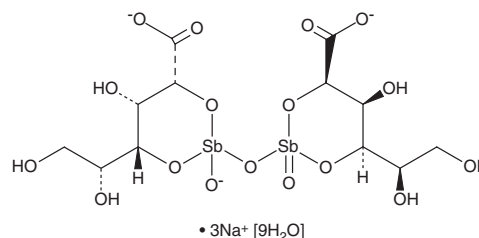
PRODUCT INFORMATION



Sodium Stibogluconate

Item No. 16038

CAS Registry No.: 16037-91-5
Formal Name: 2,4:2',4'-O-(oxydistibylidyne)bis-D-gluconic acid, trisodium salt, nonahydrate
Synonyms: Antimony Sodium Gluconate
MF: C₁₂H₁₇O₁₇Sb₂ • 3Na [9H₂O]
FW: 907.9
Purity: ≥95%
UV/Vis.: λ_{max}: 266, 321 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

Sodium stibogluconate is supplied as a crystalline solid. A stock solution may be made by dissolving the sodium stibogluconate in water. The solubility of sodium stibogluconate in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Sodium stibogluconate is a pentavalent antimony derivative that is highly active against *Leishmania* amastigotes in macrophages.¹ Sodium stibogluconate has diverse effects on both this intracellular parasite and its host cell.^{2,3}

References

1. Roberts, W.L. and Rainey, P.M. Antileishmanial activity of sodium stibogluconate fractions. *Antimicrob. Agents Chemother.* **37(9)**, 1842-1846 (1993).
2. Frézard, F., Demicheli, C., and Ribeiro, R.R. Pentavalent antimonials: New perspectives for old drugs. *Molecules* **14(7)**, 2317-2336 (2009).
3. Singh, N., Kumar, M., and Singh, R.K. Leishmaniasis: Current status of available drugs and new potential drug targets. *Asian Pac. J. Trop. Med.* **5(6)**, 485-497 (2012).

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, 10/07/2022

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