

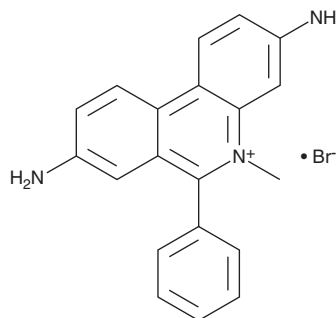
PRODUCT INFORMATION



Dimidium (bromide)

Item No. 35816

CAS Registry No.: 518-67-2
Formal Name: 3,8-diamino-5-methyl-6-phenyl-phenanthridinium, monobromide
Synonym: Phenanthridinium Compound 1553
MF: C₂₀H₁₈N₃ • Br
FW: 380.3
Purity: ≥98%
UV/Vis.: λ_{max}: 214, 295 nm
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

Dimidium (bromide) is supplied as a solid. A stock solution may be made by dissolving the dimidium (bromide) in the solvent of choice, which should be purged with an inert gas. Dimidium (bromide) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of dimidium (bromide) in these solvents is approximately 3 and 2 mg/ml, respectively.

Description

Dimidium is a trypanocide.^{1,2} *In vivo*, dimidium (0.25-0.7 mg/kg) reduces parasite burden in *T. congolense*-infected mice.¹ It also completely clears trypanosomes from the peripheral blood in zebu oxen naturally infected with *T. congolense* when administered intramuscularly at a dose of 0.75 mg/kg.² Formulations containing dimidium have previously been used in the treatment of trypanosomiasis in cattle.

References

1. Woolfe, G. The trypanocidal action of phenanthridine compounds. *Ann. Trop. Med. Parasitol.* **46(4)**, 285-288 (1952).
2. Wilde, J.K.H. Experimental use of dimidium bromide (1553) in the treatment of *T. congolense* infection in the zebu cattle. *J. Comp. Pathol. Ther.* **59(1)**, 48-53 (1949).

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, 12/13/2022

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