

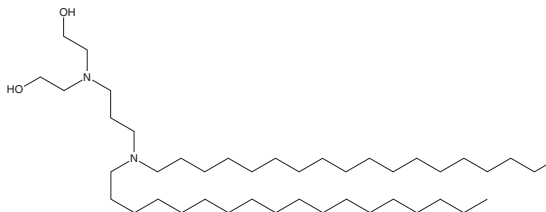
# PRODUCT INFORMATION



## Avridine

Item No. 30198

**CAS Registry No.:** 35607-20-6  
**Formal Name:** 2,2'-[[3-(dioctadecylamino)propyl]imino]bis-ethanol  
**Synonym:** CP 20,961  
**MF:** C<sub>43</sub>H<sub>90</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 667.2  
**Purity:** ≥95%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Avridine is supplied as a crystalline solid. A stock solution may be made by dissolving the avridine in the solvent of choice, which should be purged with an inert gas. Avridine is soluble in organic solvents such as ethanol and DMSO.

### Description

Avridine is a synthetic lipid amine that acts as an adjuvant.<sup>1</sup> Subplantar injection of avridine induces paw swelling, an effect that can be blocked by the non-steroidal anti-inflammatory drug (NSAID) phenylbutazone (Item No. 70400) or the alkylating agent cyclophosphamide (Item No. 13849), in rats. Avridine induces paw joint tissue destruction, bone and pannus formation, and neutrophil, fibroblast, and osteoclast infiltration in the same model. It also increases the number of CD4<sup>+</sup> T and MHC class II<sup>+</sup> cells in rat paws when administered at a dose of 1.5 mg/animal.<sup>2</sup> Avridine has been used to induce experimental rheumatoid arthritis in rats.<sup>2-4</sup>

### References

1. Chang, Y.-H., Pearson, C.M., and Abe, C. Adjuvant polyarthritis. IV. Induction by a synthetic adjuvant: Immunologic, histopathologic, and other studies. *Arthritis Rheum.* **23(1)**, 62-71 (1980).
2. Vingsbo, C., Jonsson, R., and Holmdahl, R. Avridine-induced arthritis in rats; a T cell-dependent chronic disease influenced both by MHC genes and by non-MHC genes. *Clin. Exp. Immunol.* **99(3)**, 359-363 (1995).
3. Brun, J.G., Håland, G., Haga, H.-J., *et al.* Effects of calprotectin in avridine-induced arthritis. *APMIS* **103(3)**, 233-240 (1995).
4. Holmdahl, R., Lorentzen, J.C., Lu, S., *et al.* Arthritis induced in rats with non-immunogenic adjuvants as models for rheumatoid arthritis. *Immunol. Rev.* **184**, 184-202 (2001).

#### 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, 04/10/2020

#### 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