

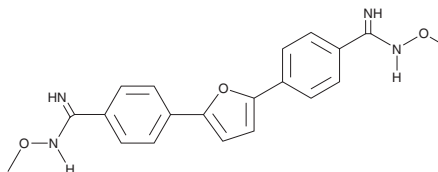
PRODUCT INFORMATION



Pafuramidine

Item No. 36669

CAS Registry No.: 186953-56-0
Formal Name: 4,4'-(2,5-furandiyl)bis[N-methoxy-benzenecarboximidamide]
Synonym: DB289
MF: C₂₀H₂₀N₄O₃
FW: 364.4
Purity: ≥90%
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

Pafuramidine is supplied as a solid. A stock solution may be made by dissolving the pafuramidine in the solvent of choice, which should be purged with an inert gas. Pafuramidine is slightly soluble in DMSO and methanol (heated).

Description

Pafuramidine is a prodrug form of the protein arginine methyltransferase 1 (PRMT1) inhibitor furamidine (Item No. 19121).¹ Oral administration of pafuramidine (25 mg/kg per day), in combination with the antiprotozoal benznidazole (Item No. 29090), reduces the number of circulating parasites and increases survival in a mouse model of *T. cruzi* infection.²

References

1. Boykin, D.W., Kumar, A., Hall, J.E., *et al.* Anti-pneumocystis activity of bis-amidoximes and bis-o-alkylamidoximes prodrugs. *Bioorganic Med. Chem. Lett.* **6(24)**, 3017-3020 (1996).
2. de Gama Jaén Batista, D., Batista, M.M., de Oliveira, G.M., *et al.* Combined treatment of heterocyclic analogues and benznidazole upon *Trypanosoma cruzi* in vivo. *PLoS One* **6(7)**, e22155 (2011).

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, 03/02/2023

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