

PRODUCT INFORMATION



Tribendimidine

Item No. 43907

CAS Registry No.: 115103-15-6

Formal Name: N',N'''-[1,4-phenylenebis(methylidynenitrilo-4,1-phenylene)]bis[N,N-dimethylethanimidamide]

MF: C₂₈H₃₂N₆

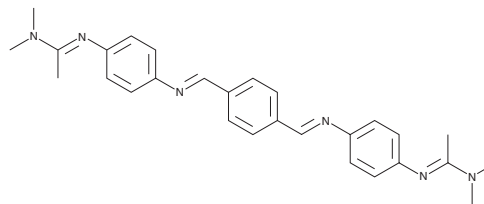
FW: 452.6

Purity: ≥98%

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

Tribendimidine is supplied as a solid. A stock solution may be made by dissolving the tribendimidine in the solvent of choice, which should be purged with an inert gas. Tribendimidine is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Tribendimidine is an anthelmintic.¹ It induces mortality in newly excysted *O. felinus* metacercariae and adult flukes (IC₅₀s = 0.23 and 0.19 μM, respectively). *In vivo*, tribendimidine (400 mg/kg) reduces worm burden in *O. felinus*-infected hamsters. Formulations containing tribendimidine have been used in the treatment of soil-transmitted helminth infections.

Reference

1. Pakharukova, M.Y., Samsonov, V.A., Serbina, E.A., et al. A study of tribendimidine effects *in vitro* and *in vivo* on the liver fluke *Opisthorchis felinus*. *Parasit. Vectors* **12**(1), 23 (2019).

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, 02/26/2026

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