

PRODUCT INFORMATION



1,3-PBIT (dihydrobromide)

Item No. 81500

CAS Registry No.: 200716-66-1

Formal Name: S,S'-1,3-phenylene-bis(1,2-ethanediyl)bis-isothiourea, dihydrobromide

MF: $C_{12}H_{18}N_4S_2 \cdot 2HBr$

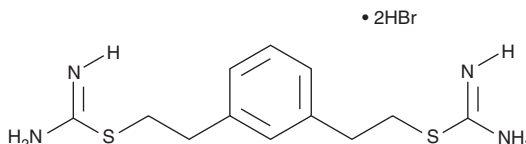
FW: 444.2

Purity: $\geq 98\%$

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Description

1,3-PBIT is a potent inhibitor of iNOS with a K_i of 47 nM compared to K_i values for eNOS and nNOS are 9 and 0.25 μM , respectively, for the purified human enzymes.¹ Its inhibition in whole cells is greatly diminished, presumably to poor membrane permeability.

Reference

1. Garvey, E.P., Oplinger, J.A., Tanoury, G.J., *et al.* Potent and selective inhibition of human nitric oxide synthases. Inhibition by non-amino acid isothioureas. *J. Biol. Chem.* **269**(43), 26669-26676 (1994).

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, 04/09/2024

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