

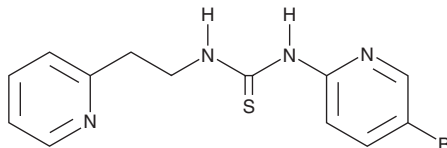
# PRODUCT INFORMATION



## Trovirdine

Item No. 43467

**CAS Registry No.:** 149488-17-5  
**Formal Name:** N-(5-bromo-2-pyridinyl)-N'-[2-(2-pyridinyl)ethyl]-thiourea  
**Synonyms:** LY300046, LY297345  
**MF:** C<sub>13</sub>H<sub>13</sub>BrN<sub>4</sub>S  
**FW:** 337.2  
**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

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

### Description

Trovirdine is a non-nucleoside reverse transcriptase inhibitor (NNRTI).<sup>1</sup> It inhibits HIV-1 wild-type reverse transcriptase (RT; IC<sub>50</sub> = 0.017 μM) and the mutants RT<sup>L100I</sup>, RT<sup>E138R</sup>, RT<sup>Y181C</sup>, RT<sup>Y188H</sup> (IC<sub>50</sub>s = 0.43, 0.22, 2.5, and 0.2 μM, respectively). Trovirdine (5 mg/kg) delays p27 antigenemia and the appearance of anti-simian immunodeficiency virus (anti-SIV) antibodies in rhesus monkeys infected with transgenic SIV expressing HIV-1 RT.<sup>2</sup> It is also active against *T. gondii* (IC<sub>50</sub> = 18.89 μM).<sup>3</sup>

### References

1. Zhang, H., Vrang, L., Bäckbro, K., *et al.* Inhibition of human immunodeficiency virus type 1 wild-type and mutant reverse transcriptases by the phenyl ethyl thiazolyl thiourea derivatives trovirdine and MSC-127. *Antiviral Res.* **28(4)**, 331-342 (1995).
2. Wang, J.-L., Elsheikha, H.M., Li, T.-T., *et al.* Efficacy of antiretroviral compounds against *Toxoplasma gondii* in vitro. *Int. J. Antimicrob. Agents* **54(6)**, 814-819 (2019).
3. Uberla, K., Stahl-Hennig, C., Böttiger, D., *et al.* Animal model for the therapy of acquired immunodeficiency syndrome with reverse transcriptase inhibitors. *Proc. Natl. Acad. Sci. USA* **92(18)**, 8210-8214 (1995).

#### 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, 05/21/2025

#### 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