

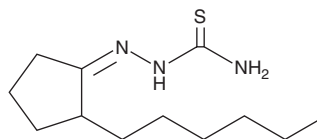
PRODUCT INFORMATION



BLT-1

Item No. 34989

CAS Registry No.: 321673-30-7
Formal Name: 2-(2-hexylcyclopentylidene)-hydrazinecarbothioamide
Synonym: Block Lipid Transport-1
MF: C₁₂H₂₃N₃S
FW: 241.4
Purity: ≥98%
UV/Vis.: λ_{max}: 273 nm
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

BLT-1 is supplied as a solid. A stock solution may be made by dissolving the BLT-1 in the solvent of choice, which should be purged with an inert gas. BLT-1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of BLT-1 in these solvents is approximately 5, 10, and 11 mg/ml, respectively.

Description

BLT-1 is an inhibitor of scavenger receptor B1 (SR-BI).¹ It inhibits uptake of [³H]cholesteryl ester-HDL in LDL receptor-deficient CHO cells expressing mouse SR-BI (IdIA[mSR-BI]) cells; IC₅₀ = 0.11 μM and the efflux of cholesterol from IdIA[mSR-BI] cells to HDL (IC₅₀ = 0.15 μM). BLT-1 (8 μM) induces developmental defects in zebrafish embryos, an effect that can be prevented by copper chloride.²

References

1. Nieland, T.J.F., Penman, M., Dori, L., *et al.* Discovery of chemical inhibitors of the selective transfer of lipids mediated by the HDL receptor SR-BI. *Proc. Natl. Acad. Sci. USA* **99(24)**, 15422-15427 (2002).
2. Raldúa, D. and Babin, P.J. BLT-1, a specific inhibitor of the HDL receptor SR-BI, induces a copper-dependent phenotype during zebrafish development. *Toxicol. Lett.* **175(1-3)**, 1-7 (2007).

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, 11/16/2022

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