

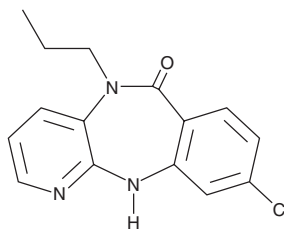
PRODUCT INFORMATION



BI-0115

Item No. 43991

CAS Registry No.: 4929-23-1
Formal Name: 9-chloro-5,11-dihydro-5-propyl-6H-pyrido[2,3-b][1,4]benzodiazepin-6-one
MF: C₁₅H₁₄ClN₃O
FW: 287.7
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

BI-0115 is supplied as a solid. A stock solution may be made by dissolving the BI-0115 in the solvent of choice, which should be purged with an inert gas. BI-0115 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

BI-0115 is an inhibitor of oxidized low-density lipoprotein receptor 1 (Ox-LDL receptor 1).¹ It inhibits oxidized LDL uptake in CHO cells expressing human Ox-LDL receptor 1 (IC₅₀ = 5.4 μM).

Reference

1. Schnapp, G., Neubauer, H., Büttner, F.H., *et al.* A small-molecule inhibitor of lectin-like oxidized LDL receptor-1 acts by stabilizing an inactive receptor tetramer state. *Commun. Chem.* **3(1)**, 75 (2020).

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, 08/19/2025

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