

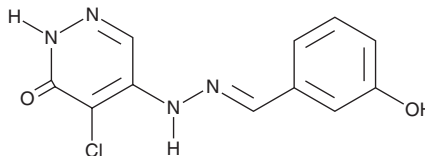
PRODUCT INFORMATION



L82-G17

Item No. 41041

CAS Registry No.: 92285-87-5
Formal Name: 4-chloro-5-(2-(3-hydroxybenzylidene)hydrazineyl)pyridazin-3(2H)-one
MF: $C_{11}H_9ClN_4O_2$
FW: 264.7
Purity: $\geq 98\%$
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

L82-G17 is supplied as a solid. A stock solution may be made by dissolving the L82-G17 in the solvent of choice, which should be purged with an inert gas. L82-G17 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

L82-G17 is an uncompetitive inhibitor of DNA ligase I.¹ It increases the retention of DNA ligase I, but not DNA ligase III, on nicked DNA-coated beads in a cell-free assay, as well as decreases the proliferation of HeLa cells when used at a concentration of 20 μM .

Reference

- Howes, T.R.L., Sallmyr, A., Brooks, R., *et al.* Structure-activity relationships among DNA ligase inhibitors: Characterization of a selective uncompetitive DNA ligase I inhibitor. *DNA Repair (Amst)* **60**, 29-39 (2017).

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, 06/03/2024

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