

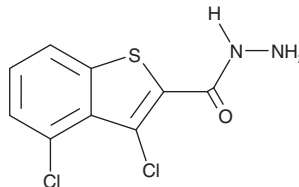
PRODUCT INFORMATION



OGG1-IN-O8

Item No. 29769

CAS Registry No.: 350997-39-6
Formal Name: 3,4-dichloro-benzo[b]thiophene-2-carboxylic acid, hydrazide
MF: C₉H₆Cl₂N₂OS
FW: 261.1
Purity: ≥98%
UV/Vis.: λ_{max}: 237, 284, 311 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

OGG1-IN-O8 is supplied as a solid. A stock solution may be made by dissolving the OGG1-IN-O8 in the solvent of choice, which should be purged with an inert gas. OGG1-IN-O8 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of OGG1-IN-O8 in these solvents is approximately 10 and 5 mg/ml, respectively.

OGG1-IN-O8 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, OGG1-IN-O8 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. OGG1-IN-O8 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

OGG1-IN-O8 is an inhibitor of 8-oxoguanine DNA glycosylase 1 (OGG1; IC₅₀ = 0.35 μM).¹ It is selective for OGG1 over Nei-like DNA glycosylase 1 (NEIL1), endonuclease III homolog 1 (NTH1), and formamidopyrimidine DNA glycosylase (Fpg) at 50 μM.

Reference

1. Lloyd, R.S., McCullough, A.K., and Donley, N. Small molecule inhibitors of 8-oxoguanine DNA glycosylase-1 (OGG1). *Oregon Health & Science University* **WO2017011834A1** (2016).

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, 10/21/2022

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