

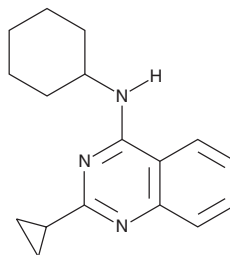
PRODUCT INFORMATION



TH10785

Item No. 43043

CAS Registry No.: 1002801-51-5
Formal Name: N-cyclohexyl-2-cyclopropyl-4-quinazolinamine
MF: C₁₇H₂₁N₃
FW: 267.4
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

TH10785 is supplied as a solid. A stock solution may be made by dissolving the TH10785 in the solvent of choice, which should be purged with an inert gas. TH10785 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

TH10785 is an 8-oxoguanine DNA glycosylase 1 (OGG1) activator.¹ It increases OGG1-mediated excision of 8-oxoguanine (8-hydroxyguanine; Item No. 89290) in duplex oligonucleotides when used at a concentration of 10 μM. TH10785 (10, 30, or 100 μM) inhibits oxidative damage in A549 cells induced by paraquat (Item No. 38521). It reduces opacity and vacuole formation induced by hydrogen peroxide (H₂O₂) in primary rat lenses when used at a concentration of 10 μM.²

References

1. Baptiste, B.A. Katchur, S.R. Fivenson, *et al.* Enhanced mitochondrial DNA repair of the common disease-associated variant, Ser326Cys, of hOGG1 through small molecule intervention. *Free Radic. Biol. Med.* **124**, 149-162 (2018)
2. Geng, W. Li, P. Zhang, G., *et al.* Targeted activation of OGG1 inhibits paraptosis in lens epithelial cells of early age-related cortical cataract. *Invest. Ophthalmol. Vis. Sci.* **66**(1), 29 (2025)

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, 04/07/2025

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