

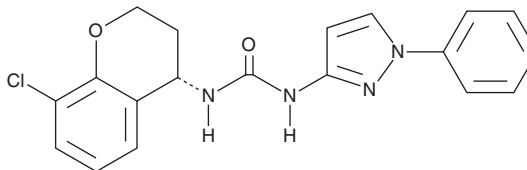
PRODUCT INFORMATION



PZL-A

Item No. 45767

CAS Registry No.: 3068830-89-4
Formal Name: N-[(4S)-8-chloro-3,4-dihydro-2H-1-benzopyran-4-yl]-N'-(1-phenyl-1H-pyrazol-3-yl)-urea
MF: C₁₉H₁₇ClN₄O₂
FW: 368.8
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

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

Description

PZL-A is an allosteric activator of mutant DNA polymerase γ (POL γ), an enzyme frequently mutated in POLG-related mitochondrial disease.¹ It stabilizes the interface between the subunits POL γ A and POL γ B in mutant POL γ and restores polymerase activity with half maximal activity concentration (AC₅₀) values of 0.16 and 0.02 μ M for POL γ ^{A467T} and POL γ ^{G848S}, respectively. PZL-A restores decreases in mitochondrial DNA (mtDNA) levels induced by ethidium bromide (Item No. 39132) and increases ATP production in fibroblasts isolated from patients with POL γ ^{A467T/G848S} or POL γ ^{W748S/R232H} mutations.

Reference

1. Valenzuela, S., Zhu, X., Macao, B., *et al.* Small molecules restore mutant mitochondrial DNA polymerase activity. *Nature* **642**(8067), 501-507 (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, 07/14/2026

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