

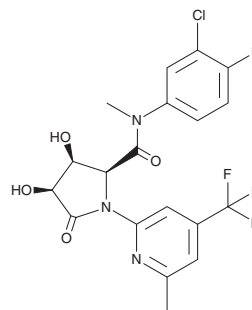
PRODUCT INFORMATION



ART812

Item No. 42904

CAS Registry No.: 2607138-82-7
Formal Name: (2S,3S,4S)-N-(3-chloro-4-fluorophenyl)-3,4-dihydroxy-N-methyl-1-[6-methyl-4-(trifluoromethyl)-2-pyridinyl]-5-oxo-2-pyrrolidinecarboxamide
MF: C₁₉H₁₆ClF₄N₃O₄
FW: 461.8
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

ART812 is supplied as a solid. A stock solution may be made by dissolving the ART812 in the solvent of choice, which should be purged with an inert gas. ART812 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

ART812 is an inhibitor of DNA polymerase θ (IC₅₀ = 7.6 nM).¹ It inhibits microhomology-mediated end joining (MMEJ) in a reporter assay using HEK293 cells (EC₅₀ = 240 nM). ART812 (150 mg/kg twice per day) increases blood levels of micronucleated reticulocytes in mice. It inhibits tumor development in a *Brca1*- and *Shld2*-knockout MDA-MB-436 breast cancer rat xenograft model of tumor initiation when administered at a dose of 100 mg/kg per day.

Reference

1. Stockley, M.L., Benedetti, G., Blencowe, P., *et al.* Discovery, characterization, and structure-based optimization of small-molecule in vitro and in vivo probes for human DNA polymerase theta. *J. Med. Chem.* **65**(20), 13879-13891 (2022).

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, 02/19/2025

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