

PRODUCT INFORMATION



TP-021

Item No. 41315

CAS Registry No.: 2130878-25-8

Formal Name: 6-[[2-chloro-4-nitro-5-[[tetrahydro-2H-pyran-4-yl]oxy]phenyl]amino]-3,4-dihydro-2(1H)-quinolinone

MF: C₂₀H₂₀ClN₃O₅

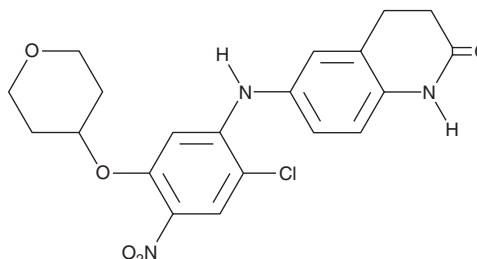
FW: 417.8

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

TP-021 is supplied as a solid. A stock solution may be made by dissolving the TP-021 in the solvent of choice, which should be purged with an inert gas. TP-021 is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

Description

TP-021 is an inhibitor of the protein-protein interaction between Bcl-6 and Bcl-6 corepressor (BCoR).¹ It inhibits the protein-protein interaction between the Bcl-6 BTB domain and BCoR in a mammalian two-hybrid assay (IC₅₀ = 0.72 μM).

Reference

1. Yasui, T., Yamamoto, T., Sakai, N., *et al.* Discovery of a novel B-cell lymphoma 6 (BCL6)-corepressor interaction inhibitor by utilizing structure-based drug design. *Bioorg. Med. Chem.* **25**(17), 4876-4886 (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, 07/01/2024

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