

PRODUCT INFORMATION

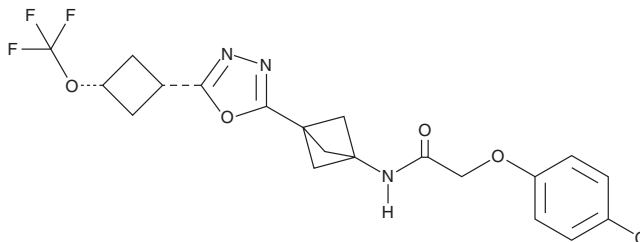


DNL343

Item No. 41674

CAS Registry No.: 2278265-85-1
Formal Name: 2-(4-chlorophenoxy)-N-[3-[5-[*cis*-3-(trifluoromethoxy)cyclobutyl]-1,3,4-oxadiazol-2-yl]bicyclo[1.1.1]pent-1-yl]-acetamide

MF: C₂₀H₁₉ClF₃N₃O₄
FW: 457.8
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

DNL343 is supplied as a solid. A stock solution may be made by dissolving the DNL343 in the solvent of choice, which should be purged with an inert gas. DNL343 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

DNL343 is an activator of eukaryotic translation initiation factor 2B (eIF2B; IC₅₀ = 9.8 nM in a reporter assay) and a derivative of 2BAct (Item No. 37788).¹ It inhibits sodium arsenite-induced formation of stress granules positive for TAR DNA-binding protein 43 (TDP-43) in an *in vitro* model of amyotrophic lateral sclerosis (ALS) using H4 cells (IC₅₀ = 12.46 nM). DNL343 (50 mg/kg) decreases the expression of integrated stress response (ISR) genes in Eif2b5^{R191H} mutant mice in a model of neurodegeneration.

Reference

1. Craig, R.A.I., De Vicente, J., Estrada, A.A., *et al.* Discovery of DNL343: A potent, selective, and brain-penetrant eIF2B activator designed for the treatment of neurodegenerative diseases. *J. Med. Chem.* **67**(7), 5758-5782 (2024).

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, 09/17/2024

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