

PRODUCT INFORMATION

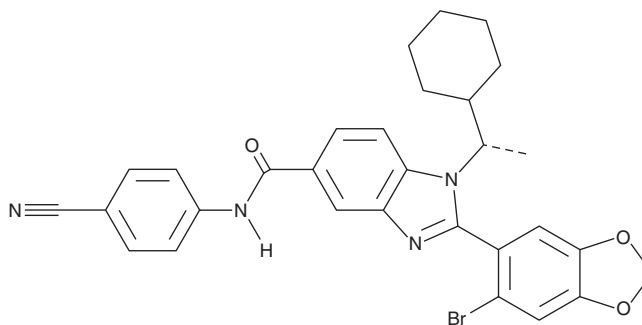


AZ3451

Item No. 29671

CAS Registry No.: 2100284-59-9
Formal Name: 2-(6-bromo-1,3-benzodioxol-5-yl)-N-(4-cyanophenyl)-1-[(1S)-1-cyclohexylethyl]-1H-benzimidazole-5-carboxamide

MF: C₃₀H₂₇BrN₄O₃
FW: 571.5
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 298 nm
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

AZ3451 is supplied as a solid. A stock solution may be made by dissolving the AZ3451 in the solvent of choice, which should be purged with an inert gas. AZ3451 is soluble in DMSO.

Description

AZ3451 is an allosteric antagonist of proteinase-activated receptor 2 (PAR2) that inhibits calcium mobilization by wild-type and 10 mutant forms of PAR2 induced by the PAR2 ligand SLIGRL in a FLIPR assay (IC₅₀s = 23 and 12-780 nM, respectively).¹ It is selective for PAR2 over PAR4 and PAR1 (IC₅₀s = <2.5, 380, and >50,000 nM, respectively, in a β-arrestin-2 recruitment assay). AZ3451 completely inhibits SLIGRL-induced phosphorylation of ERK in 1321N1 astrocytoma cells when used at a concentration of 10 μM.

Reference

1. Cheng, R.K.Y., Fiez-Vandal, C., Schlenker, O., *et al.* Structural insight into allosteric modulation of protease-activated receptor 2. *Nature* **545**(7652), 112-115 (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, 11/01/2022

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