

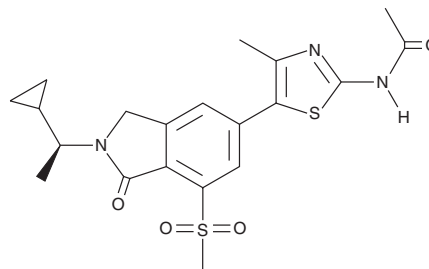
PRODUCT INFORMATION



AZD 3458

Item No. 42104

CAS Registry No.: 2132961-46-5
Formal Name: N-[5-[2-[(1S)-1-cyclopropylethyl]-2,3-dihydro-7-(methylsulfonyl)-1-oxo-1H-isoindol-5-yl]-4-methyl-2-thiazolyl]-acetamide
MF: C₂₀H₂₃N₃O₄S₂
FW: 433.5
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

AZD 3458 is supplied as a solid. A stock solution may be made by dissolving the AZD 3458 in the solvent of choice, which should be purged with an inert gas. AZD 3458 is sparingly soluble (1-10 mg/ml) in methanol and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

AZD 3458 is an inhibitor of PI3K γ (IC₅₀ = 0.79 nM).¹ It is selective for PI3K γ over PI3K α , PI3K β , and PI3K δ (IC₅₀s = 7.94, >31.62, and 0.316 μ M, respectively), PI3K δ 2 α , PI3K δ 2 γ , and PI3K δ 3 (IC₅₀s = >10, 3.16, and 7.94 μ M, respectively), and a panel of 395 kinases at 1 μ M but does not inhibit PI3K δ 2 β (IC₅₀ = 31.62 nM). AZD 3458 (0.1, 0.5, and 2.5 mg/kg) reduces LPS-induced increases in the number of neutrophils in bronchoalveolar lavage fluid (BALF) in a rat model of acute lung inflammation.

Reference

1. Pemberton, N., Mogemark, M., Arlbrandt, S., *et al.* Discovery of highly isoform selective orally bioavailable phosphoinositide 3-kinase (PI3K)- γ inhibitors. *J. Med. Chem.* **61**(12), 5435-5441 (2018).

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, 11/05/2024

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