

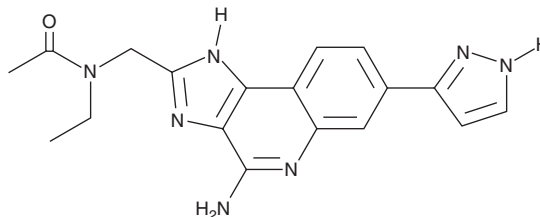
PRODUCT INFORMATION



BMS 986299

Item No. 40652

CAS Registry No.: 2242952-69-6
Formal Name: N-[[4-amino-7-(1H-pyrazol-3-yl)-3H-imidazo[4,5-c]quinolin-2-yl]methyl]-N-ethyl-acetamide
MF: C₁₈H₁₉N₇O
FW: 349.4
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

BMS 986299 is supplied as a solid. A stock solution may be made by dissolving the BMS 986299 in the solvent of choice, which should be purged with an inert gas. BMS 986299 is slightly soluble (0.1-1 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of BMS 986299 can be prepared by directly dissolving the solid in aqueous buffers. BMS 986299 is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

BMS 986299 is an agonist of the NOD-like receptor protein 3 (NLRP3) inflammasome.¹ It induces IL-1 β secretion from mouse macrophages expressing the human NLRP3 inflammasome (EC₅₀ = 1.28 μ M).

Reference

1. Glick, G., Ghosh, S., Roush, W.R., *et al.* Substituted imidazo-quinolines as NLRP3 modulators. *Innate Tumor Immunity, Inc.* **WO2018/152396A1** (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, 05/01/2024

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