

PRODUCT INFORMATION

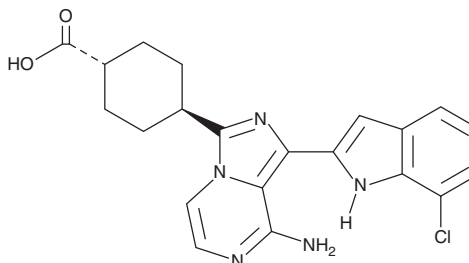


OXA-01

Item No. 20456

CAS Registry No.: 936889-68-8
Formal Name: *trans*-4-[8-amino-1-(7-chloro-1H-indol-2-yl)imidazo[1,5-a]pyrazin-3-yl]-cyclohexanecarboxylic acid

MF: C₂₁H₂₀ClN₅O₂
FW: 409.9
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 336 nm
Supplied as: A crystalline 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

OXA-01 is supplied as a crystalline solid. A stock solution may be made by dissolving the OXA-01 in the solvent of choice. OXA-01 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of OXA-01 in these solvents is approximately 3 and 1 mg/ml, respectively.

OXA-01 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, OXA-01 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. OXA-01 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

OXA-01 is an inhibitor of both mTORC1 and mTORC2 (IC₅₀s = 11 and 29 nM, respectively).¹ It is selective for mTOR over PI3K isoforms and DNA protein kinase.¹ OXA-01 reduces VEGF production in tumors in association with decreased vessel sprouting.¹

Reference

1. Falcon, B.L., Barr, S., Gokhale, P.C., *et al.* Reduced VEGF production, angiogenesis, and vascular regrowth contribute to the antitumor properties of dual mTORC1/mTORC2 inhibitors. *Cancer Res.* **71(5)**, 1573-1583 (2011).

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, 10/20/2022

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