

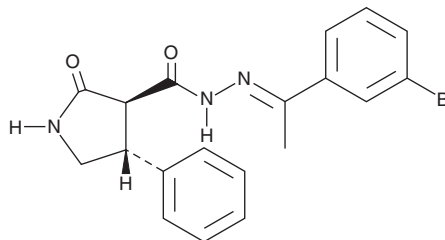
PRODUCT INFORMATION



AC-264613

Item No. 28777

CAS Registry No.: 1051487-82-1
Formal Name: (3R,4S)-rel-2-oxo-4-phenyl-3-pyrrolidinecarboxylic acid, 2-[1-(3-bromophenyl)ethylidene]hydrazide
MF: C₁₉H₁₈BrN₃O₂
FW: 400.3
Purity: ≥95%
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

AC-264613 is supplied as a solid. A stock solution may be made by dissolving the AC-264613 in the solvent of choice, which should be purged with an inert gas. AC-264613 is soluble in the organic solvent DMSO at a concentration of approximately 40 mg/ml.

Description

AC-264613 is an agonist of proteinase-activated receptor 2 (PAR2).¹ It induces calcium mobilization, phosphatidylinositol hydrolysis, and cell proliferation in NIH3T3 cells expressing PAR2 (EC₅₀s = 100, 125.8, and 31.6 nM, respectively). AC-264613 (3 µg/paw) induces paw edema and thermal hyperalgesia in rats.

Reference

1. Gardell, L.R., Ma, J.N., Seitzberg, J.G., *et al.* Identification and characterization of novel small-molecule protease-activated receptor 2 agonists. *J. Pharmacol. Exp. Ther.* **327**(3), 799-808 (2008).

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, 11/08/2022

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