

PRODUCT INFORMATION

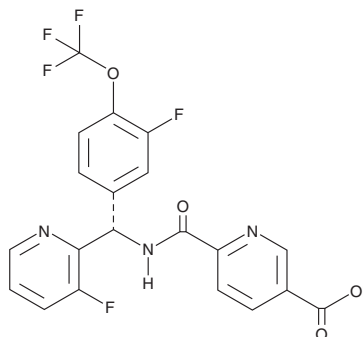


AMG 333

Item No. 36825

CAS Registry No.: 1416799-28-4
Formal Name: 6-[[[(S)-(3-fluoro-2-pyridinyl)[3-fluoro-4-(trifluoromethoxy)phenyl]methyl]amino]carbonyl]-3-pyridinecarboxylic acid

MF: C₂₀H₁₂F₅N₃O₄
FW: 453.3
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 268 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

AMG 333 is supplied as a solid. A stock solution may be made by dissolving the AMG 333 in the solvent of choice, which should be purged with an inert gas. AMG 333 is soluble in DMSO.

Description

AMG 333 is an antagonist of transient receptor potential melastatin 8 (TRPM8; IC₅₀s = 13 and 20 for the human and rat channels, respectively).¹ It is selective for TRPM8 over TRP vanilloid 1 (TRPV1), TRPV3, TRPV4, and TRP ankyrin 1 (TRPA1; IC₅₀s = >20, >20, >20, and >40 μM, respectively), as well as a panel of 100 kinases and 144 additional targets at 1 and 10 μM, respectively. AMG 333 (0.6-3 mg/kg) decreases wet-dog shakes, a behavioral marker of cold sensation, induced by the TRPM8 agonist icilin (Item No. 10137) in rats.

Reference

1. Horne, D.B., Biswas, K., Brown, J., *et al.* Discovery of TRPM8 antagonist (S)-6-(((3-fluoro-4-(trifluoromethoxy)phenyl)(3-fluoropyridin-2-yl)methyl)carbamoyl)nicotinic acid (AMG 333), a clinical candidate for the treatment of migraine. *J. Med. Chem.* **61**(18), 8186-8201 (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, 11/30/2022

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