

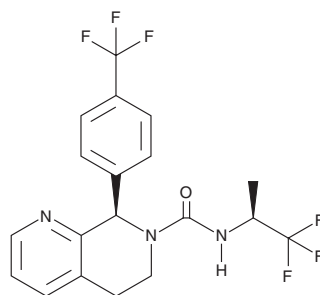
PRODUCT INFORMATION



AMG 2850

Item No. 42906

CAS Registry No.: 1470018-52-0
Formal Name: 5,8R-dihydro-N-[(1S)-2,2,2-trifluoro-1-methylethyl]-8-[4-(trifluoromethyl)phenyl]-1,7-naphthyridine-7(6H)-carboxamide
MF: C₁₉H₁₇F₆N₃O
FW: 417.4
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

AMG 2850 is supplied as a solid. A stock solution may be made by dissolving the AMG 2850 in the solvent of choice, which should be purged with an inert gas. AMG 2850 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

AMG 2850 is an antagonist of transient receptor potential melastatin 8 (TRPM8).¹ It inhibits menthol- or cold-induced intracellular calcium release in CHO cells expressing rat TRPM8 (IC₅₀s = 7.3 and 150 nM, respectively). AMG 2850 is selective for TRPM8 over transient receptor potential ankyrin 1 (TRPA1), transient receptor potential vanilloid 1 (TRPV1), TRPV3, and TRPV4 (IC₅₀s = >20, >10, >10, and >10 μM, respectively). *In vivo*, AMG 2850 (100 mg/kg) decreases body temperature in rats and mice. It also inhibits wet-dog shakes induced by icilin (Item No. 10137) in rats when administered at a dose of 10 mg/kg.²

References

1. Gavva, N.R., Davis, C., Lehto, S.G., *et al.* Transient receptor potential melastatin 8 (TRPM8) channels are involved in body temperature regulation. *Mol. Pain.* **8**, 36 (2012).
2. Lehto, S.G., Weyer, A.D., Zhang, M., *et al.* AMG2850, a potent and selective TRPM8 antagonist, is not effective in rat models of inflammatory mechanical hypersensitivity and neuropathic tactile allodynia. *Naunyn Schmiedebergs Arch. Pharmacol.* **388**(4), 465-476 (2015).

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, 03/06/2025

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