

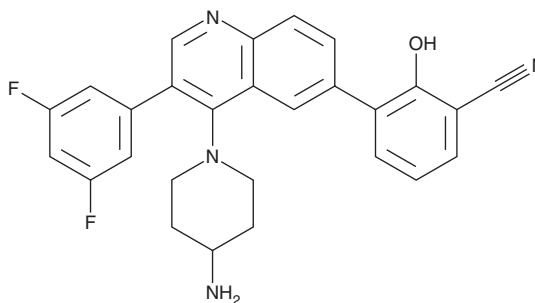
PRODUCT INFORMATION



Paltusotine

Item No. 42041

CAS Registry No.: 2172870-89-0
Formal Name: 3-[4-(4-amino-1-piperidiny)-3-(3,5-difluorophenyl)-6-quinolinyl]-2-hydroxy-benzonitrile
Synonym: CRN00808
MF: C₂₇H₂₂F₂N₄O
FW: 456.5
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

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

Description

Paltusotine is a nonpeptide somatostatin receptor 2 (SST₂) agonist.¹ It selectively inhibits cAMP production in cells expressing SST₂ over cells expressing SST₁, SST₃, SST₄, or SST₅ (EC₅₀s = 25, >1,000, >1,000, >1,000, and >1,000 nM, respectively). *In vivo*, paltusotine (3, 10, and 30 mg/kg) reduces growth hormone release stimulated by growth hormone-releasing hormone (GHRH) in rats. It also reduces plasma levels of insulin-like growth factor 1 (IGF-1) in beagle dogs.

Reference

1. Zhao, J., Wang, S., Markison, S., *et al.* Discovery of paltusotine (CRN00808), a potent, selective, and orally bioavailable non-peptide SST₂ agonist. *ACS Med. Chem. Lett.* **14**(1), 66-74 (2022).

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, 11/11/2024

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