

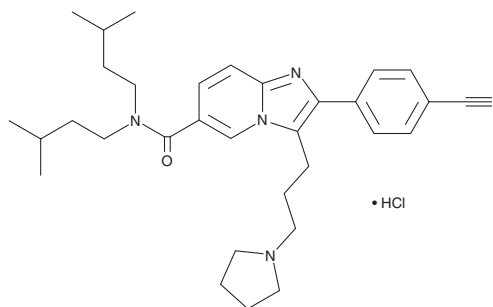
PRODUCT INFORMATION



SNT-207707 (hydrochloride)

Item No. 41524

CAS Registry No.: 1064662-40-3
Formal Name: 2-(4-cyanophenyl)-N,N-bis(3-methylbutyl)-3-[3-(1-pyrrolidinyl)propyl]-imidazo[1,2-a]pyridine-6-carboxamide, monohydrochloride
MF: C₃₂H₄₃N₅O • HCl
FW: 550.2
Purity: ≥98%
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

SNT-207707 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the SNT-207707 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SNT-207707 (hydrochloride) is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

SNT-207707 is a melanocortin receptor 4 (MC4R) antagonist (IC₅₀ = 5 nM).¹ It increases food intake in mice when administered at a dose of 20 mg/kg. SNT-207707 (30 mg/kg per day) inhibits tumor-induced weight loss without inhibiting tumor growth in a C26 murine colon carcinoma model of tumor-induced cachexia.

Reference

1. Weyermann, P., Dallmann, R., Magyar, J., *et al.* Orally available selective melanocortin-4 receptor antagonists stimulate food intake and reduce cancer-induced cachexia in mice. *PLoS One* **4**(3), e4774 (2009).

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, 08/08/2024

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