

PRODUCT INFORMATION



MK-3207

Item No. 42490

CAS Registry No.: 957118-49-9

Formal Name: 8R-(3,5-difluorophenyl)-10-oxo-N-[(2R)-1,1',2',3-tetrahydro-2'-oxospiro[2H-indene-2,3'-[3H]pyrrolo[2,3-b]pyridin]-5-yl]-6,9-diazaspiro[4.5]decane-9-acetamide

MF: C₃₁H₂₉F₂N₅O₃

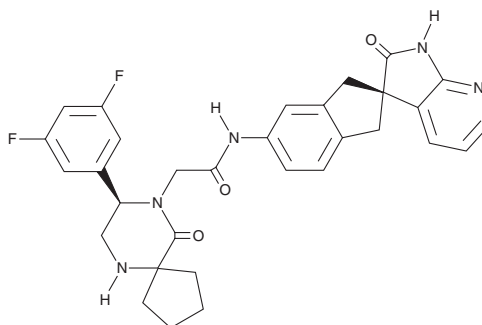
FW: 557.6

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

MK-3207 is supplied as a solid. A stock solution may be made by dissolving the MK-3207 in the solvent of choice, which should be purged with an inert gas. MK-3207 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

MK-3207 is an antagonist of the calcitonin gene-related peptide (CGRP) receptor ($K_i = 22$ pM for the recombinant human receptor).¹ It is selective for the CGRP receptor over adrenomedullin receptor 1 (AM₁) and AM₂, calcitonin (CT) receptor, and amylin 3 receptor (AMY₃; K_i s = 16.5, 0.156, 1.9, and 0.128 μ M, respectively) but does inhibit AMY₁ to a lesser extent ($K_i = 750$ pM). MK-3207 inhibits CGRP-induced cAMP production in HEK293 cells expressing the human CGRP receptor ($IC_{50} = 120$ pM). It decreases capsaicin-induced forearm dermal vasodilation in rhesus monkeys in a dose-dependent manner.

Reference

1. Salvatore, C.A., Moore, E.L., Calamari, A., et al. Pharmacological properties of MK-3207, a potent and orally active calcitonin gene-related peptide receptor antagonist. *J. Pharmacol. Exp. Ther.* **333**(1), 152-160 (2010).

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, 12/02/2025

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