

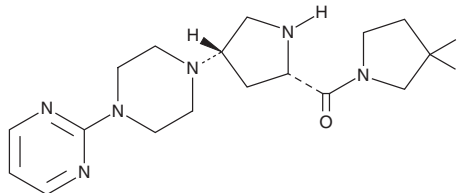
PRODUCT INFORMATION



Gosogliptin

Item No. 33220

CAS Registry No.: 869490-23-3
Formal Name: (3,3-difluoro-1-pyrrolidinyl)[(2S,4S)-4-[4-(2-pyrimidinyl)-1-piperazinyl]-2-pyrrolidinyl]-methanone
Synonym: PF 734200
MF: C₁₇H₂₄F₂N₆O
FW: 366.4
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

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

Description

Gosogliptin is an inhibitor of dipeptidyl peptidase 4 (DPP-4; IC₅₀ = 0.013 μM).¹ It is selective for DPP-4 over DPP-2, -3, -8, and -9 (IC₅₀s = 3.3, >30, 7, and 5.98 μM, respectively), the serine proteases APP, FAP, and POP (IC₅₀s = >30, 10.3, and >30 μM, respectively), and a panel of enzymes, receptors, and ion channels (IC₅₀s = >10 μM for all).

Reference

1. Ammirati, M.J., Andrews, K.M., Boyer, D.D., *et al.* (3,3-Difluoro-pyrrolidin-1-yl)-[(2S,4S)-(4-(4-pyrimidin-2-yl-piperazin-1-yl)-pyrrolidin-2-yl)]-methanone: A potent, selective, orally active dipeptidyl peptidase IV inhibitor. *Bioorg. Med. Chem. Lett.* **19(7)**, 1991-1995 (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, 12/09/2022

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