

PRODUCT INFORMATION



5-hydroxy Saxagliptin (hydrochloride)

Item No. 26710

Formal Name: (1S,3S,5S)-2-((S)-2-amino-2-((1r,3R,5S,7S)-3,5-dihydroxyadamantan-1-yl)acetyl)-2-azabicyclo[3.1.0]hexane-3-carbonitrile, monohydrochloride

MF: C₁₈H₂₅N₃O₃ • HCl

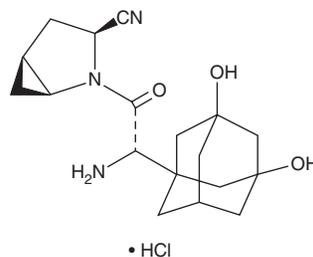
FW: 367.9

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

5-hydroxy Saxagliptin (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the 5-hydroxy saxagliptin (hydrochloride) in the solvent of choice. 5-hydroxy Saxagliptin (hydrochloride) is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

5-hydroxy Saxagliptin is a major active metabolite of saxagliptin (Item No. 23697).¹ 5-hydroxy Saxagliptin inhibits dipeptidyl peptidase 4 (DPP-4; K_i = 2.6 nM for human recombinant enzyme). It is selective for DPP-4 over DPP-8 and DPP-9 (K_{i,s} = 2,495 and 423 nM, respectively, for human recombinant enzymes).

Reference

1. Wang, A., Dorso, C., Kopcho, L., *et al.* Potency, selectivity and prolonged binding of saxagliptin to DPP4: Maintenance of DPP4 inhibition by saxagliptin in vitro and ex vivo when compared to a rapidly-dissociating DPP4 inhibitor. *BMC Pharmacol.* **12**, 2 (2012).

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, 11/03/2022

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