

Sitagliptin

Catalog #: 27106

Lot #: 111121

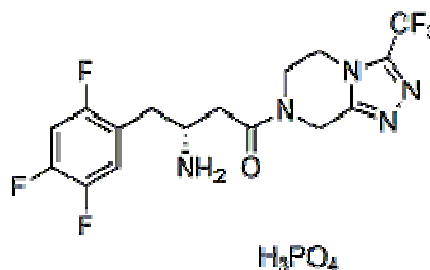
Size: 200 mg

Structure:

CAS Registry #: 654671-78-0

Purity: >99%

Chemical Formula: C₁₆H₁₅F₆N₅O.H₃PO₄



Molecular Weight: 505.31

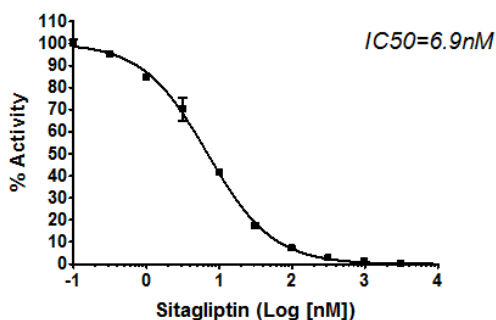
Description: Sitagliptin, also known as MK-0431 and Januvia, is a Dipeptidyl Peptidase IV Inhibitor.

Appearance: White to beige powder.

Solubility: Soluble in DMSO.

Biological Activity: Sitagliptin as a DPP-4 inhibitor has IC₅₀ value of 6.9 nM when testing treatment of patients with type 2 diabetes mellitus.

DPP4 Activity



Storage/Stability: Store at or below -20 °C.

Quality Control: The inhibitor was determined by NMR analysis.

Reference(s):

1. Gallwitz B. Drugs Today (Barc). 2007 Nov; **43(11)**:801-14
2. Su-Jin Kim et al. Diabetes, 2010 July; **59(7)**:1739-1750

For R&D use only. Not for human, drug, diagnostic, household or other uses. MSDS available. Product of USA.
6044 Cornerstone Court West, Suite E, San Diego, CA 92121

Tel: (858) 829-3082 • Fax: (858) 481-8694

Email: info@bpsbioscience.com • Visit: www.bpsbioscience.com • © 2011 BPS Bioscience Inc.