

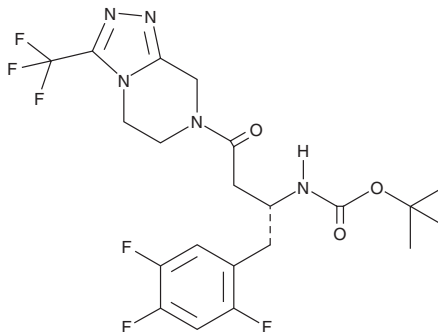
PRODUCT INFORMATION



N-Boc Sitagliptin

Item No. 35312

CAS Registry No.: 486460-23-5
Formal Name: N-[(1R)-3-[5,6-dihydro-3-(trifluoromethyl)-1,2,4-triazolo[4,3-a]pyrazin-7(8H)-yl]-3-oxo-1-[(2,4,5-trifluorophenyl)methyl]propyl]-carbamic acid, 1,1-dimethylethyl ester
MF: C₂₁H₂₃F₆N₅O₃
FW: 507.4
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

N-Boc sitagliptin is supplied as a solid. A stock solution may be made by dissolving the N-boc sitagliptin in the solvent of choice, which should be purged with an inert gas. N-Boc sitagliptin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N-boc sitagliptin in DMSO is approximately 16 mg/ml and approximately 20 mg/ml in DMF. N-Boc sitagliptin is slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of N-boc sitagliptin can be prepared by directly dissolving the solid in aqueous buffers. The solubility of N-boc sitagliptin in PBS (pH 7.2) is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N-Boc sitagliptin is an intermediate in the synthesis of the dipeptidyl peptidase 4 (DPP-4) inhibitor (-)-sitagliptin (Item No. 13252).¹

Reference

1. Anjibabu, R., Boggu, J.M.R., Shekhar, P., *et al.* Stereoselective synthesis of dipeptidyl peptidase-4 (DPP-4) inhibitor, (R)-sitagliptin. *ChemistrySelect* **1**(17), 5445-5447 (2016).

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, 12/12/2022

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