

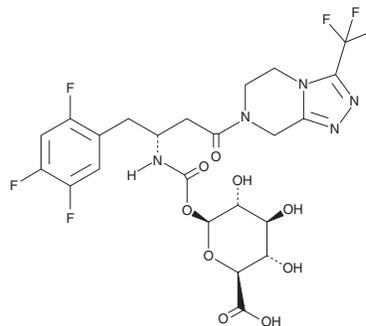
PRODUCT INFORMATION



(-)-Sitagliptin Carbamoyl Glucuronide

Item No. 26629

CAS Registry No.: 940002-59-5
Formal Name: 1-[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] carbamate] β-D-glucopyranuronic acid
Synonym: (R)-Sitagliptin Carbamoyl Glucuronide
MF: C₂₃H₂₃F₆N₅O₉
FW: 627.5
Purity: ≥90%
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

(-)-Sitagliptin carbamoyl glucuronide is supplied as a solid. A stock solution may be made by dissolving the (-)-sitagliptin carbamoyl glucuronide in water. We do not recommend storing the aqueous solution for more than one day.

Description

(-)-Sitagliptin carbamoyl glucuronide is a minor phase II metabolite of the dipeptidyl peptidase 4 (DPP-4) inhibitor (-)-sitagliptin (Item No. 13252).¹ (-)-Sitagliptin carbamoyl glucuronide is formed by N-carbamoyl glucuronidation and has been found in rat and dog plasma following administration of (-)-sitagliptin.

Reference

1. Beconi, M.G., Reed, J.R., Teffera, Y., *et al.* Disposition of the dipeptidyl peptidase 4 inhibitor sitagliptin in rats and dogs. *Drug Metab. Dispos.* **35**(4), 525-532 (2007).

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, 12/21/2022

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