

PRODUCT INFORMATION

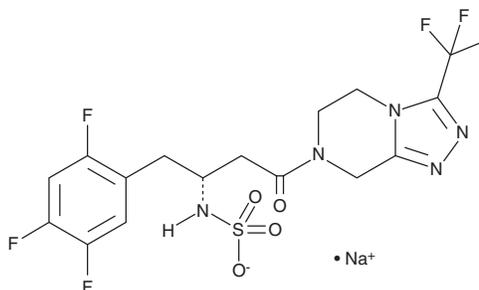


Sitagliptin N-Sulfate (sodium salt)

Item No. 26625

CAS Registry No.: 2470126-25-9
Formal Name: (R)-(4-oxo-4-(3-(trifluoromethyl)-5,6-dihydro-[1,2,4]triazolo[4,3-a]pyrazin-7(8H)-yl)-1-(2,4,5-trifluorophenyl)butan-2-yl)sulfamate, monosodium salt

MF: C₁₆H₁₄F₆N₅O₄S • Na
FW: 509.4
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

Sitagliptin N-sulfate (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the sitagliptin N-sulfate (sodium salt) in the solvent of choice, which should be purged with an inert gas. Sitagliptin N-sulfate (sodium salt) is slightly soluble in methanol.

Description

Sitagliptin N-sulfate is a phase II metabolite of the dipeptidyl peptidase 4 (DPP-4) inhibitor (-)-sitagliptin (Item No. 13252).¹ Sitagliptin N-sulfate is formed *via* sulfation and has been found in rat bile following oral 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, 01/27/2023

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