

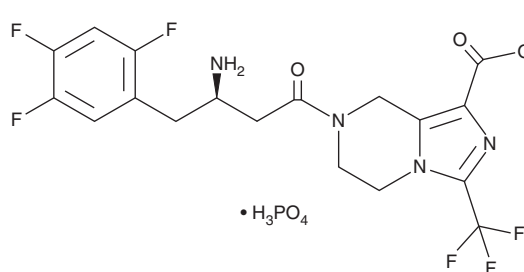
PRODUCT INFORMATION



Retagliptin (phosphate)

Item No. 45052

CAS Registry No.: 1256756-88-3
Formal Name: 7-[(3R)-3-amino-1-oxo-4-(2,4,5-trifluorophenyl)butyl]-5,6,7,8-tetrahydro-3-(trifluoromethyl)-2 imidazo[1,5-a]pyrazine-1-carboxylic acid, methyl ester, monophosphate
Synonym: SP2086
MF: $C_{19}H_{18}F_6N_4O_3 \cdot H_3PO_4$
FW: 562.4
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Retagliptin (phosphate) is supplied as a solid. A stock solution may be made by dissolving the retagliptin (phosphate) in the solvent of choice, which should be purged with an inert gas. Retagliptin (phosphate) slightly soluble (0.1-1 mg/ml) in DMSO.

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 retagliptin (phosphate) can be prepared by directly dissolving the solid in aqueous buffers. Retagliptin (phosphate) is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Retagliptin is an inhibitor of dipeptidyl peptidase 4 (DPP-4).¹ It induces autophagy and inhibits melanogenesis induced by α -melanocyte-stimulating hormone (α -MSH) in B16-F1 melanocytes when used at concentrations of 10 and 100 μM .

Reference

1. Kim, S.H., Bae, J.-E., Park, N., *et al.* Evaluation of teneligliptin and retagliptin on the clearance of melanosome by melanophagy in B16F1 cells. *Cosmetics* **11**(2), 35 (2024).

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, 02/06/2026

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