

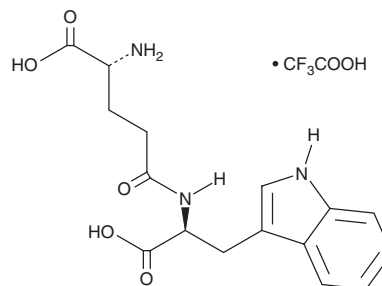
PRODUCT INFORMATION



Golotimod (trifluoroacetate salt)

Item No. 40833

CAS Registry No.: 2828433-07-2
Formal Name: D-γ-glutamyl-L-tryptophan, 2,2,2-trifluoroacetate
Synonyms: γ-D-Glutamyl-L-Tryptophan, SCV-07
MF: C₁₆H₁₉N₃O₅ • CF₃COOH
FW: 447.4
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

Golotimod (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the golotimod (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Golotimod (trifluoroacetate salt) is soluble (≥ 10 mg/ml) in ethanol and 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 golotimod (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Golotimod (trifluoroacetate salt) is soluble in (≥ 10 mg/ml) PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Golotimod is a synthetic dipeptide.¹ *In vivo*, golotimod (100 µg/kg per day) decreases the incidence of vaginal infection by herpes simplex virus 2 (HSV-2) and cumulative mean lesion scores in HSV-2-infected guinea pigs.

Reference

1. Rose II, W.A., Tuthill, C., and Pyles, R.B. An immunomodulating dipeptide, SCV-07, is a potential therapeutic for recurrent genital herpes simplex virus type 2 (HSV-2). *Int. J. Antimicrob. Agents* **32**(3), 262-266 (2008).

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, 06/20/2024

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