

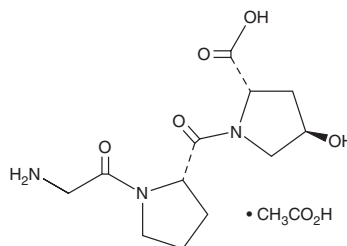
PRODUCT INFORMATION



H-Gly-Pro-Hyp-OH (acetate)

Item No. 35390

CAS Registry No.: 103192-50-3
Formal Name: glycyl-L-prolyl-4R-hydroxy-L-proline, monoacetate
Synonyms: H-GP-Hyp-OH, Tripeptide-29
MF: $C_{12}H_{19}N_3O_5 \cdot C_2H_4O_2$
FW: 345.3
Purity: $\geq 95\%$
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

H-Gly-Pro-Hyp-OH (acetate) is supplied as a solid. Aqueous solutions of H-Gly-Pro-Hyp-OH (acetate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of H-Gly-Pro-Hyp-OH (acetate) in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

H-Gly-Pro-Hyp-OH is a peptide dipeptidyl peptidase 4 (DPP-4) inhibitor ($IC_{50} = 2.51 \text{ mM}$).¹

Reference

1. Hatanaka, T., Kawakami, K., and Uraji, M. Inhibitory effect of collagen-derived tripeptides on dipeptidylpeptidase-IV activity. *J. Enzyme Inhib. Med. Chem.* **29**(6), 823-828 (2014).

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