

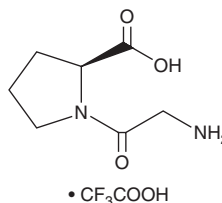
PRODUCT INFORMATION



Glycyl-L-Proline (trifluoroacetate salt)

Item No. 36182

CAS Registry No.: 76397-13-2
Synonym: Glycylproline
MF: $C_7H_{12}N_2O_3 \cdot CF_3COOH$
FW: 286.2
Purity: $\geq 95\%$
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

Glycyl-L-proline (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the glycyl-L-proline (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Glycyl-L-proline (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of glycyl-L-proline (trifluoroacetate salt) in these solvents is approximately 1, 5, and 2 mg/ml, respectively.

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 glycyl-L-proline (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of glycyl-L-proline (trifluoroacetate salt) in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Glycyl-L-proline is a dipeptide composed of glycine and L-proline (Item No. 30772).¹ It has been used in the study of dipeptide transport in enterocyte basolateral membrane and brush-membrane border vesicles.^{1,2}

References

1. Rajendran, V.M., Harig, J.M., and Ramaswamy, K. Characteristics of glycyl-L-proline transport in intestinal brush-border membrane vesicles. *Am. J. Physiol.* **252**(2 Pt. 1), G281-6 (1987).
2. Dyer, J., Beechey, R.B., Gorvel, J.P., et al. Glycyl-L-proline transport in rabbit enterocyte basolateral-membrane vesicles. *Biochem. J.* **269**(3), 565-571 (1990).

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, 12/19/2022

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