

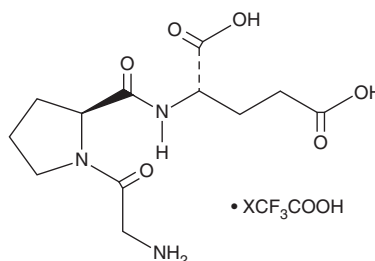
# PRODUCT INFORMATION



## Gly-Pro-Glu (trifluoroacetate salt)

Item No. 43926

**Formal Name:** glycyl-L-prolyl-L-glutamic acid, trifluoroacetate salt  
**Synonyms:** Gly-Pro-Glu-OH, GPE, IGF-1 (1-3), Insulin-like Growth Factor 1 (1-3)  
**Peptide Sequence:** GPE-OH  
**MF:** C<sub>12</sub>H<sub>19</sub>N<sub>3</sub>O<sub>6</sub> • XCF<sub>3</sub>COOH  
**FW:** 301.3  
**Purity:** ≥98%  
**Supplied as:** A low-melting 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

Gly-Pro-Glu (trifluoroacetate salt) is supplied as a low-melting solid. A stock solution may be made by dissolving the Gly-Pro-Glu (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Gly-Pro-Glu (trifluoroacetate salt) is sparingly soluble (1-10 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 Gly-Pro-Glu (trifluoroacetate salt) can be prepared by directly dissolving the low-melting solid in aqueous buffers. Gly-Pro-Glu (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Gly-Pro-Glu is a peptide antagonist of the NMDA receptor (IC<sub>50</sub> = 170 μM).<sup>1</sup> It potentiates potassium-induced release of acetylcholine (ACh; Item No. 23829) in rat cortical slices and dopamine (Item Nos. 36532 | 21992) in rat striatal slices when used at concentrations of 0.1 and 10 nM. Gly-Pro-Glu (100 μM) increases the proliferation of primary human Müller glial cells.<sup>2</sup> It increases the proliferation of primary mouse embryonic neural stem cells and induces migration of the same cells in a wound healing assay when used at a concentration of 100 μM.<sup>3</sup> Gly-Pro-Glu increases locomotor activity in the open field test in aged rats when administered during the neonatal period at a dose of 30 μg/animal for 12 days.<sup>4</sup>

### References

1. Sara, V.R., Carlsson-Skwirut, C., Bergman, T., *et al.* Identification of Gly-Pro-Glu (GPE), the aminoterminal tripeptide of insulin-like growth factor 1 which is truncated in brain, as a novel neuroactive peptide. *Biochem. Biophys. Res. Commun.* **165**(2), 766-771 (1989).
2. Ikeda, T., Waldbillig, R.J., and Puro, D.G. Truncation of IGF-I yields two mitogens for retinal Müller glial cells. *Brain Res.* **686**(1), 87-92 (1995).
3. Almengló, C., Devesa, P., Devesa, J., *et al.* GPE promotes the proliferation and migration of mouse embryonic neural stem cells and their progeny in vitro. *Int. J. Mol. Sci.* **18**(6), 1280 (2017).
4. Sara, V.R., Carlsson-Skwirut, C., Drakenberg, K., *et al.* The biological role of truncated insulin-like growth factor-1 and the tripeptide GPE in the central nervous system. *Ann. N.Y. Acad. Sci.* **692**, 183-191 (1993).

#### 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, 08/11/2025

#### 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