

# PRODUCT INFORMATION



## GHRP-2 (trifluoroacetate salt)

Item No. 28811

**Formal Name:** D-alanyl-3-(2-naphthalenyl)-D-alanyl-L-alanyl-L-tryptophyl-D-phenylalanyl-L-lysineamide, 2,2,2-trifluoroacetate

**Synonyms:** D-Ala-3-(2-naphthalenyl)-D-Ala-L-Ala-L-Trp-D-Phe-L-lysineamide, Growth Hormone-Releasing Peptide 2, KP-102, Pralmorelin

**MF:** C<sub>45</sub>H<sub>55</sub>N<sub>9</sub>O<sub>6</sub> • CF<sub>3</sub>COOH

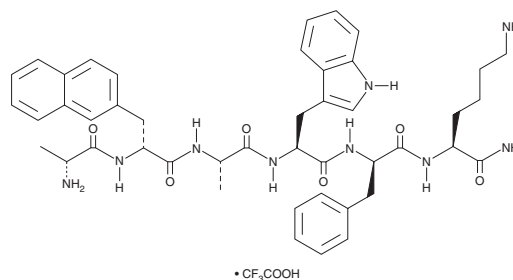
**FW:** 932.0

**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

GHRP-2 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the GHRP-2 (trifluoroacetate salt) in water. The solubility of GHRP-2 (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

GHRP-2 is an orally bioavailable synthetic growth hormone (GH) secretagogue and an agonist of the GH secretagogue receptor (GHS-R), which is also known as the ghrelin receptor.<sup>1</sup> It induces GH release from primary rat pituitary cells (EC<sub>50</sub> = 1.8 nM) and binds to human hypothalamic membranes (IC<sub>50</sub> = 16 nM).<sup>2,3</sup> GHRP-2 increases blood GH levels in anesthetized rats (ED<sub>50</sub> = 30 nmol/kg) and serum GH levels in conscious pigs (ED<sub>50</sub> = 0.6 nmol/kg).<sup>2</sup> It also increases food intake and body weight gain in rats when administered at a dose of 100 µg/kg per day.<sup>4</sup>

### References

1. Peroni, C.N., Hayashida, C.Y., Nascimento, N., *et al.* Growth hormone response to growth hormone-releasing peptide-2 in growth hormone-deficient little mice. *Clinics (Sao Paulo)* **67(3)**, 265-272 (2012).
2. Ankersen, M., Johansen, N.L., Madsen, K., *et al.* A new series of highly potent growth hormone-releasing peptides derived from ipamorelin. *J. Med. Chem.* **41(19)**, 3699-3704 (1998).
3. Muccioli, G., Ghè, C., Ghigo, M.C., *et al.* Specific receptors for synthetic GH secretagogues in the human brain and pituitary gland. *J. Endocrinol.* **157(1)**, 99-106 (1998).
4. Granado, M., Priego, T., Martín, A.I., *et al.* Anti-inflammatory effect of the ghrelin agonist growth hormone-releasing peptide-2 (GHRP-2) in arthritic rats. *Am. J. Physiol. Endocrinol. Metab.* **288(3)**, E486-E492 (2005).

#### 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, 12/05/2022

#### 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