

# PRODUCT INFORMATION



## Peptide YY (24-36) (trifluoroacetate salt)

Item No. 45087

**Formal Name:** L-leucyl-L-arginyl-L-histidyl-L-tyrosyl-L-leucyl-L-asparaginyl-L-leucyl-L-valyl-L-threonyl-L-arginyl-L-glutamyl-L-arginyl-L-tyrosinamide, trifluoroacetate salt

**Synonyms:** Leu-Arg-His-Tyr-Leu-Asn-Leu-Val-Thr-Arg-Gln-Arg-Tyr-NH<sub>2</sub>, PYY (24-36)

**Peptide Sequence:** LRHYLNLVTRQRY-NH<sub>2</sub>

**MF:** C<sub>78</sub>H<sub>127</sub>N<sub>27</sub>O<sub>18</sub> • XCF<sub>3</sub>COOH

**FW:** 1,731.0

**Purity:** ≥98%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

H—Leu—Arg—His—Tyr—Leu—Asn—Leu—Val—Thr—Arg—Gln—Arg—Tyr—NH<sub>2</sub>  
• XCF<sub>3</sub>COOH

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Peptide YY (24-36) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the peptide YY (24-36) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Peptide YY (24-36) (trifluoroacetate salt) is soluble (≥10 mg/ml) in 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 peptide YY (24-36) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Peptide YY (24-36) (trifluoroacetate salt) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Peptide YY (24-36) is a synthetic C-terminal peptide fragment of the anorectic peptide hormone peptide YY (Item No. 24776).<sup>1</sup> Unlike peptide YY, peptide YY (24-36) does not inhibit cholecystokinin-8-induced pancreatic protein secretion in dogs.

### Reference

1. Yoshinaga, K., Mochizuki, T., Yanaihara, N., *et al.* Structural requirements of peptide YY for biological activity at enteric sites. *Am. J. Physiol.* **263**(5 Pt 1), G695-G701 (1992).

**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, 02/25/2026

### 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