

PRODUCT INFORMATION



Peptide YY (human) (trifluoroacetate salt)

Item No. 24776

Formal Name: L-tyrosyl-L-prolyl-L-isoleucyl-L-lysyl-L-prolyl-L- α -glutamyl-L-alanyl-L-prolylglycyl-L- α -glutamyl-L- α -aspartyl-L-alanyl-L-seryl-L-prolyl-L- α -glutamyl-L- α -glutamyl-L-leucyl-L-asparaginy-L-arginyl-L-tyrosyl-L-tyrosyl-L-alanyl-L-seryl-L-leucyl-L-arginyl-L-histidyl-L-tyrosyl-L-leucyl-L-asparaginy-L-leucyl-L-valyl-L-threonyl-L-arginyl-L-glutaminy-L-arginyl-L-tyrosinamide, trifluoroacetate salt

Tyr-Pro-Ile-Lys-Pro-Glu-Ala-Pro-Gly-Glu-
Asp-Ala-Ser-Pro-Glu-Glu-Leu-Asn-Arg-Tyr-
Tyr-Ala-Ser-Leu-Arg-His-Tyr-Leu-Asn-Leu-
Val-Thr-Arg-Gln-Arg-Tyr-NH₂

• XCF₃COOH

Synonyms: Peptide Tyrosine Tyrosine, PYY

MF: C₁₉₄H₂₉₅N₅₅O₅₇ • XCF₃COOH

FW: 4,309.8

Purity: ≥95%

Supplied as: A lyophilized powder

Storage: -20°C

Stability: ≥4 years

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

Laboratory Procedures

Peptide YY (PYY) (human) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the PYY (human) (trifluoroacetate salt) in water. The solubility of PYY (human) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PYY is a 36-amino acid peptide and anorectic gut hormone agonist for the neuropeptide Y receptors Y₁, Y₂, Y₅, and Y₆ with EC₅₀ values of 0.7, 0.58, 1, and 0.8 nM, respectively, for suppression of forskolin-induced cAMP accumulation.^{1,2} PYY is cleaved *in vivo* to PYY (3-36) (Item No. 16502). Release of PYY occurs postprandially in proportion to calorie intake from intestinal enteroendocrine L cells, indicating it may be a satiety signal.³ In humans, PYY inhibits food intake and gastrointestinal transit in both lean and obese subjects.

References

1. Gerald, C., Walker, M.W., Criscione, L., *et al.* A receptor subtype involved in neuropeptide-Y-induced food intake. *Nature* **382(6587)**, 168-171 (1996).
2. Mullins, D.E., Guzzi, M., Xia, L., *et al.* Pharmacological characterization of the cloned neuropeptide Y y₆ receptor. *Eur. J. Pharmacol.* **395(2)**, 87-93 (2000).
3. Holzer, P., Reichmann, F., and Farzi, A. Neuropeptide Y, peptide YY and pancreatic polypeptide in the gut-brain axis. *Neuropeptides* **46(6)**, 261-274 (2012).

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