

# PRODUCT INFORMATION



## PACAP-related Peptide (rat) (trifluoroacetate salt)

Item No. 24773

CAS Registry No.: 132769-35-8

**Formal Name:** L- $\alpha$ -aspartyl-L-valyl-L-alanyl-L-histidyl-L- $\alpha$ -glutamyl-L-isoleucyl-L-leucyl-L-asparaginyl-L- $\alpha$ -glutamyl-L-alanyl-L-tyrosyl-L-arginyl-L-lysyl-L-valyl-L-leucyl-L- $\alpha$ -aspartyl-L-glutamyl-L-leucyl-L-seryl-L-alanyl-L-arginyl-L-lysyl-L-tyrosyl-L-leucyl-L-glutamyl-L-seryl-L-methionyl-L-valyl-L-alanine, trifluoroacetate salt

H — Asp — Val — Ala — His — Glu — Ile — Leu — Asn — Glu — Ala —

Tyr — Arg — Lys — Val — Leu — Asp — Gln — Leu — Ser — Ala —

Arg — Lys — Tyr — Leu — Gln — Ser — Met — Val — Ala — NH<sub>2</sub>

• XCF<sub>3</sub>COOH

**Synonyms:** Pituitary Adenylate Cyclase-activating Peptide-related Peptide (rat), PRP (rat)

**MF:** C<sub>148</sub>H<sub>242</sub>N<sub>42</sub>O<sub>45</sub>S • XCF<sub>3</sub>COOH

**FW:** 3,361.8

**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

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

### Description

PRP is an endogenous 29-amino acid peptide that belongs to the secretin/glucagon superfamily of peptides, which includes secretin (Item No. 24561), glucagon (Item No. 24204), glucagon-like peptide-1 (GLP-1; Item No. 24460), GLP-2 (Item No. 24414), and pituitary adenylate cyclase-activating polypeptide (PACAP; Item Nos. 24769 | 24770).<sup>1,2</sup> It is expressed in rat hypothalamus as well as within the nerves of the median eminence, the anterior pituitary, bed nucleus of the stria terminalis, cerebellum, cerebral cortex, and amygdala.<sup>3,4</sup> PRP is also expressed in vaginal, uterine cervical, uterine horn, fallopian tube, and ovarian tissues from the rat female genital tract and is present in extracts of male testis tissue.<sup>5,6</sup>

### References

1. Vaudry, D., Falluel-Morel, A., Bourgault, S., et al. *Pharmacol. Rev.* **61**(3), 283-357 (2009).
2. Tam, J.K., Lee, L.T., and Chow, B.K. *Peptides* **28**(9), 1920-1929 (2007).
3. Hannibal, J., Mikkelsen, J.D., Clausen, H., et al. *Regul. Pept.* **55**(2), 133-148 (1995).
4. Mikkelsen, J.D., Hannibal, J., Fahrenkrug, J., et al. *J. Neuroendocrinol.* **7**(1), 47-55 (1995).
5. Fahrenkrug, J. and Hannibal, J. *Neuroscience* **73**(4), 1049-1060 (1996).
6. Hannibal, J. and Fahrenkrug, J. *Regul. Pept.* **55**(1), 111-115 (1995).

#### 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, 11/18/2022

#### 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](http://WWW.CAYMANCHEM.COM)