

# PRODUCT INFORMATION



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

Item No. 24772

**Formal Name:** L- $\alpha$ -aspartyl-L-valyl-L-alanyl-L-histidylglycyl-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-lysyl-L-histidyl-L-leucyl-L-glutamyl-L-seryl-L-leucyl-L-valyl-L-alanine, trifluoroacetate salt

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

**MF:**  $C_{139}H_{229}N_{41}O_{42} \cdot XCF_3COOH$

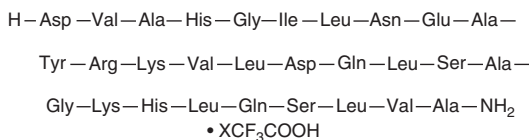
**FW:** 3,146.6

**Purity:**  $\geq 95\%$

**Supplied as:** A lyophilized powder

**Storage:**  $-20^{\circ}C$

**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

PACAP-related peptide (PRP) (human) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the PRP (human) (trifluoroacetate salt) in water. The solubility of PRP (human) (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 normal human pancreas and adrenal gland tissue and in some tumors that produce vasoactive intestinal peptide (VIP).<sup>3</sup> PRP (1-29) is secreted by CHO-K1 cells that express human recombinant preproPACAP.<sup>4</sup>

### References

1. Vaudry, D., Falluel-Morel, A., Bourgault, S., *et al.* Pituitary adenylate cyclase-activating polypeptide and its receptors: 20 years after the discovery. *Pharmacol. Rev.* **61**(3), 283-357 (2009).
2. Tam, J.K., Lee, L.T., and Chow, B.K. PACAP-related peptide (PRP)--molecular evolution and potential functions. *Peptides* **28**(9), 1920-1929 (2007).
3. Fahrenkrug, J., Buhl, T., and Hannibal, J. PreproPACAP-derived peptides occur in VIP-producing tumours and co-exist with VIP. *Regul. Pept.* **58**(3), 89-98 (1995).
4. Okazaki, K., Kimura, C., Kosaka, T., *et al.* Expression of human pituitary adenylate cyclase activating polypeptide (PACAP) cDNA in CHO cells and characterization of the products. *FEBS Lett.* **298**(1), 49-56 (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, 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)