

PRODUCT INFORMATION



Acetyl PACAP (1-38) (human, mouse, ovine, porcine, rat) (trifluoroacetate salt) Item No. 24774

Formal Name: N-acetyl-L-histidyl-L-seryl-L- α -aspartylglycyl-L-isoleucyl-L-phenylalanyl-L-threonyl-L- α -aspartyl-L-seryl-L-tyrosyl-L-seryl-L-arginyl-L-tyrosyl-L-arginyl-L-lysyl-L-glutamyl-L-methionyl-L-alanyl-L-valyl-L-lysyl-L-lysyl-L-tyrosyl-L-leucyl-L-alanyl-L-alanyl-L-valyl-L-leucylglycyl-L-lysyl-L-arginyl-L-tyrosyl-L-lysyl-L-glutamyl-L-arginyl-L-valyl-L-lysyl-L-asparaginyl-L-lysineamide, trifluoroacetate salt



Synonym: Acetyl Pituitary Adenylate Cyclase-activating Peptide (1-38)

MF: $\text{C}_{205}\text{H}_{333}\text{N}_{63}\text{O}_{54}\text{S} \cdot \text{XCF}_3\text{COOH}$

• XCF_3COOH

FW: 4,576.3

Purity: $\geq 95\%$

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

Acetyl PACAP (1-38) (human, mouse, ovine, porcine, rat) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the acetyl PACAP (1-38) (human, mouse, ovine, porcine, rat) (trifluoroacetate salt) in water. The solubility of acetyl PACAP (1-38) (human, mouse, ovine, porcine, rat) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Acetyl pituitary adenylate cyclase-activating peptide (PACAP) (1-38) is an N-terminally acetylated form of PACAP (1-38) (Item No. 24770).¹ It binds to the PACAP receptor PAC_1 ($\text{IC}_{50} = 5.6$ nM for the human receptor) and induces calcium mobilization and proliferation of PC12 cells ($\text{EC}_{50}\text{s} = 4.4$ and 0.83 nM, respectively) with an efficacy comparable to PACAP (1-38). Acetyl PACAP (1-38) is resistant to degradation by dipeptidyl peptidase 4 (DPP-4) with a 10-fold longer half-life than PACAP (1-38), respectively.

Reference

1. Bourgault, S., Vaudry, D., Botia, B., et al. Novel stable PACAP analogs with potent activity towards the PAC1 receptor. *Peptides* **29(6)**, 919-932 (2008).

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