

PRODUCT INFORMATION

PEN (human) (acetate)

Item No. 43151

Formal Name: L-alanyl-L-alanyl-L- α -aspartyl-L-histidyl-L- α -aspartyl-L-valylglycyl-L-seryl-L- α -glutamyl-L-leucyl-L-prolyl-L-prolyl-L- α -glutamylglycyl-L-valyl-L-leucylglycyl-L-alanyl-L-leucyl-L-leucyl-L-arginyl-L-valine, acetate

Peptide Sequence: AADHDVGSELPPEGVLGALLRV-OH

MF: $C_{97}H_{159}N_{27}O_{32} \cdot XC_2H_4O_2$

FW: 2,215.5

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years

H—Ala—Ala—Asp—His—Asp—Val—Gly—Ser—Glu—Leu—
Pro—Pro—Glu—Gly—Val—Leu—Gly—Ala—Leu—Leu—
Arg—Val—OH

$\cdot XCH_3CO_2H$

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

Laboratory Procedures

PEN (human) (acetate) is supplied as a solid. Aqueous solutions of PEN (human) (acetate) can be prepared by directly dissolving the solid in aqueous buffers. PEN (human) (acetate) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

PEN is a neuropeptide and a fragment of ProSAAS, a pro-protein that is cleaved into smaller peptides that regulate neurohormone activity.¹ PEN was originally identified as a ligand for GPR83, a G protein-coupled receptor (GPCR) involved in regulating food consumption. However, it does not induce calcium mobilization in HEK293 cells expressing human GPR83.

Reference

1. Giesecke, Y., Asimi, V., Stulberg, V., et al. Is the neuropeptide PEN a ligand of GPR83? *Int. J. Mol. Sci.* **24**(20), 15117 (2023).

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, 05/08/2025

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