

PRODUCT INFORMATION



BigLEN (mouse) (acetate)

Item No. 36766

Formal Name: L-leucyl-L- α -glutamyl-L-asparaginyl-L-prolyl-L-seryl-L-prolyl-L-glutamyl-L-alanyl-L-prolyl-L-alanyl-L-arginyl-L-arginyl-L-leucyl-L-leucyl-L-prolyl-L-proline, acetate

Peptide Sequence: LENPSPQAPARRLLPP-OH

MF: $C_{78}H_{130}N_{24}O_{22} \cdot XC_2H_4O_2$

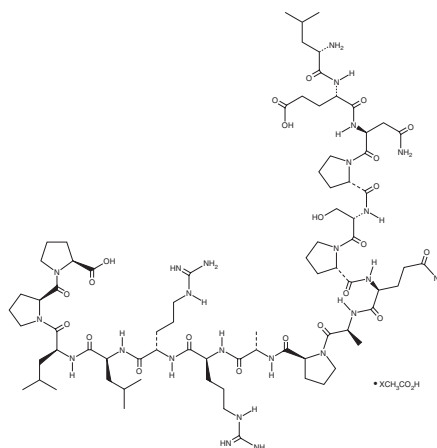
FW: 1,756.0

Purity: $\geq 90\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years



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

Laboratory Procedures

BigLEN (mouse) (acetate) is supplied as a solid. A stock solution may be made by dissolving the BigLEN (mouse) (acetate) in water. We do not recommend storing the aqueous solution for more than one day.

Description

BigLEN is a neuropeptide and fragment of the pro-protein ProSAAS, involved in food intake, and an agonist of the orphan G protein-coupled receptor 171 (GPR171).¹ It selectively increases intracellular calcium levels in CHO cells expressing GPR171 over those expressing GPR19, GPR108, or GPR165 at 1 μM . BigLEN (10 μM) increases neurite outgrowth in mouse Neuro2a neuroblastoma cells.

Reference

1. Gomes, I., Aryal, D.K., Wardman, J.H., *et al.* GPR171 is a hypothalamic G protein-coupled receptor for BigLEN, a neuropeptide involved in feeding. *Proc. Natl. Acad. Sci. USA* **110**(40), 16211-16216 (2013).

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, 08/17/2023

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