Enterostatin (human)

Reduces food intake

This pentapeptide, APGPR, has a number of physiological activities, including inhibition of insulin secretion, suppression of fat intake, and memory enhancement. *In vivo*, it is generated by the cleavage of its precursor protein, procolipase.

Ordering Information

Order Online »

BML-P234-0001 1mg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Lyophilized solid.

CAS 117830-79-2

497 MW

Purity ≥95%

Sequence Ala-Pro-Gly-Pro-Arg

Solubility Soluble in Water.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com