

PRODUCT INFORMATION



BRINP2-related Peptide

Item No. 43361

Overview and Properties

Contents:	This vial contains 1 or 5 mg of lyophilized peptide trifluoroacetate salt.	
Formal Name:	L-threonyl-L-histidyl-L-arginyl-L-isoleucyl-L-leucyl-L-arginyl-L-arginyl-L-leucyl-L-phenylalanyl-L-asparaginyl-L-leucyl-L-cysteinamide	
Synonym:	BRP	H—Thr—His—Arg—Ile—Leu—Arg—Arg—Leu—Phe—Asn—
Peptide Sequence:	THRILRRLFNLC	Leu—Cys—NH ₂
MF:	C ₆₈ H ₁₁₇ N ₂₅ O ₁₄ S	
FW:	1,540.9	
Form:	Lyophilized	
Storage:	-20°C (as supplied)	
Stability:	≥1 year	
Storage Buffer:	Soluble in water or phosphate buffer up to 30 mg/ml	

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

Description

BRINP2-related peptide is an anorexigenic peptide.¹ It is formed from BRINP2 via cleavage by proconvertases and is expressed in the brain and cerebrospinal fluid (CSF). *In vivo*, BRINP2-related peptide (5 mg/kg) reduces food intake and increases the time between meals in mice. It reduces fasting glucose and insulin levels, subcutaneous white adipose tissue (WAT), brown adipose tissue (BAT), liver mass, and body weight without reducing lean mass in a mouse model of high-fat diet-induced obesity and diabetes, independent of the glucagon-like peptide 1 receptor (GLP-1R), leptin, or melanocortin receptor 4 (MC4R). Intracerebroventricular administration of BRINP2-related peptide (0.5 mg/kg) reduces food intake and increases *Fos* expression in the dorsomedial hypothalamic nucleus, preoptic hypothalamic nucleus, tuberal nucleus, and arcuate nucleus, but not the hindbrain or brainstem, in mice.

Reference

1. Coassolo, L., Danneskiold-Samsøe, N.B., Nguyen, Q., *et al.* Prohormone cleavage prediction uncovers a non-incretin anti-obesity peptide. *Nature* (2025).

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, 07/15/2026

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