

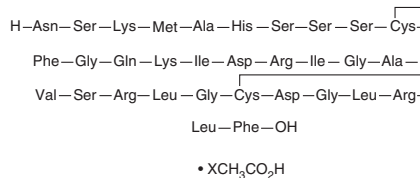
PRODUCT INFORMATION



Brain Natriuretic Peptide-32 (rat) (acetate)

Item No. 39724

Synonyms: B-type Natriuretic Peptide-32, BNP-32, BNP (14-45), Brain Natriuretic Polypeptide-32
Peptide Sequence: NSKMAHSSSCFGQKIDRIGAVSRLGCDGLRLF-OH
MF: C₁₄₆H₂₃₉N₄₇O₄₄S₃ • XC₂H₄O₂
FW: 3,452.9
Purity: ≥98%
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

Brain Natriuretic Peptide-32 (BNP-32) (rat) (acetate) is supplied as a solid. A stock solution may be made by dissolving the BNP-32 (rat) (acetate) in water. We do not recommend storing the aqueous solution for more than one day.

Description

BNP-32 is a peptide fragment of the cardiovascular hormone precursor proBNP.¹ It binds to natriuretic peptide receptors on rat vascular smooth muscle cells (IC₅₀ = 7.3 nM) and induces cGMP accumulation in the same cells (EC₅₀ = 170 nM). BNP-32 induces relaxation of norepinephrine-precontracted isolated rat aortic strips (EC₅₀ = 6.7 nM). It increases urine and urinary electrolyte excretion and induces hypotension in rats when administered at a dose of 3 µg/kg.

Reference

1. Furuya, M., Ohnuma, N., Takehisa, M., *et al.* Pharmacological activities of brain natriuretic peptides of human, porcine and rat origin. *Eur. J. Pharmacol.* **200(2-3)**, 233-237 (1991).

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, 10/19/2023

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