

PRODUCT INFORMATION

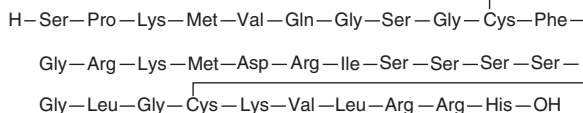


Nesiritide (acetate)

Item No. 34532

CAS Registry No.: 1684439-46-0

Formal Name: cyclic (10→26)-disulfide, L-seryl-L-prolyl-L-lysyl-L-methionyl-L-valyl-L-glutaminyglycyl-L-serylglycyl-L-cysteinyl-L-phenylalanylglycyl-L-arginyl-L-lysyl-L-methionyl-L-α-aspartyl-L-arginyl-L-isoleucyl-L-seryl-L-seryl-L-seryl-L-serylglycyl-L-leucylglycyl-L-cysteinyl-L-lysyl-L-valyl-L-leucyl-L-arginyl-L-arginyl-L-histidine, monoacetate



• CH₃CO₂H

MF: C₁₄₃H₂₄₄N₅₀O₄₂S₄ • C₂H₄O₂

FW: 3,524.1

Purity: ≥95%

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

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

Description

Nesiritide is a 32-amino acid natriuretic peptide and recombinant form of human brain natriuretic peptide (Item No. 24541).¹ It increases *ex vivo* vasorelaxation induced by the calcium ionophore A23187 (Item No. 22030) in porcine hearts isolated from a model of acute coronary occlusion when administered as a 2 µg/kg bolus dose followed by 0.01 µg/kg per minute infusion.² Formulations containing nesiritide have previously been used in the treatment of congestive heart failure.

References

1. Hobbs, R.E. and Mills, R.M. Therapeutic potential of nesiritide (recombinant b-type natriuretic peptide) in the treatment of heart failure. *Expert Opin. Investig. Drugs* **8**(7), 1063-1072 (1999).
2. Lazar, H.L., Bao, Y., Siwik, D., *et al.* Nesiritide enhances myocardial protection during the revascularization of acutely ischemic myocardium. *J. Card. Surg.* **24**(5), 600-605 (2009).

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, 11/22/2022

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