

# PRODUCT INFORMATION



## AOD-9604 (acetate)

Item No. 36790

**Formal Name:** cyclic (7→14)-disulfide,  
L-tyrosyl-L-leucyl-L-arginyl-L-  
isoleucyl-L-valyl-L-glutamyl-  
L-cysteinyl-L-arginyl-L-seryl-L-  
valyl-L-α-glutamylglycyl-L-seryl-  
L-cysteinylglycyl-L-phenylalanine,  
acetate

**MF:** C<sub>78</sub>H<sub>123</sub>N<sub>23</sub>O<sub>23</sub>S<sub>2</sub> • XC<sub>2</sub>H<sub>4</sub>O<sub>2</sub>

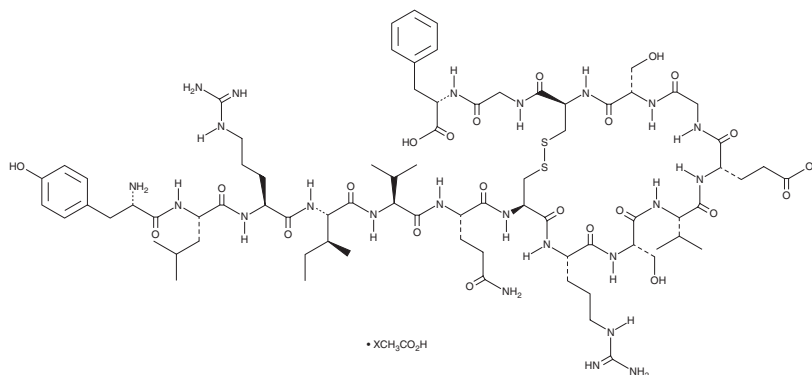
**FW:** 1,815.1

**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

AOD-9604 (acetate) is supplied as a solid. A stock solution may be made by dissolving the AOD-9604 (acetate) in the solvent of choice, which should be purged with an inert gas. AOD-9604 (acetate) is slightly soluble in DMSO.

### Description

AOD-9604 is a synthetic lipolytic 15-amino acid peptide fragment derived from human growth hormone.<sup>1,2</sup> *In vivo*, AOD-9604 (250 µg/kg per day) decreases body weight, epididymal adipose tissue mass, and brown adipose tissue weight in *ob/ob* mice.<sup>1</sup> It also enhances cartilage regeneration in a rabbit model of collagenase-induced knee osteoarthritis.<sup>2</sup>

### References

- Heffernan, M., Summers, R.J., Thorburn, A., *et al.* The effects of human GH and its lipolytic fragment (AOD9604) on lipid metabolism following chronic treatment in obese mice and β<sub>3</sub>-AR knock-out mice. *Endocrinology* **142**(12), 182-189 (2001).
- Kwon, D.R. and Park, G.Y. Effect of intra-articular injection of AOD9604 with or without hyaluronic acid in rabbit osteoarthritis model. *Ann. Clin. Lab. Sci.* **45**(4), 426-432 (2015).

#### 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, 12/06/2022

#### 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