

PRODUCT INFORMATION



Oligopeptide-20 (acetate)

Item No. 45612

Formal Name: L-arginyl-L-arginyl-L-leucyl-L- α -glutamyl-L-methionyl-L-tyrosyl-L-cysteinyl-L-alanyl-L-prolyl-L-leucyl-L-lysyl-L-proline

Peptide Sequence: RRLEMYCAPLKP-OH

MF: $C_{65}H_{109}N_{19}O_{16}S_2 \cdot XC_2H_4O_2$

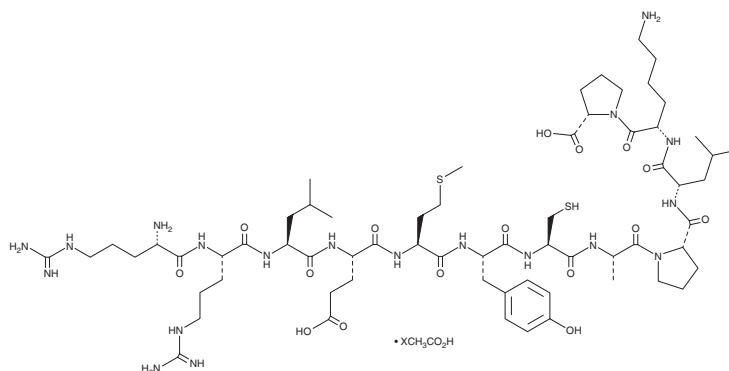
FW: 1,476.8

Purity: $\geq 95\%$

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

Oligopeptide-20 (acetate) is supplied as a solid. Aqueous solutions of oligopeptide-20 (acetate) can be prepared by directly dissolving the solid in aqueous buffers. Oligopeptide-20 (acetate) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Oligopeptide-20 is a peptide that corresponds to amino acids 55-66 of mature insulin-like growth factor 1 (IGF-1).¹ It increases growth of keratinocytes and fibroblasts and the levels of procollagen, laminin, and fibronectin in keratinocytes. Oligopeptide-20 (100 mg/animal), as part of a peptide nanosome-based skin cream formulation, increases thickness of the keratinocyte and epidermal skin layer in mice.

Reference

1. Chung, Y.J., Kim, Y.D., and Kim, E.M. Peptides having activities of insulin like growth factor-1 and their uses. **WO2007/102686A1** (2007).

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, 06/02/2026

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