

PRODUCT INFORMATION



Oligopeptide-41 (acetate)

Item No. 45628

Formal Name: (4S,7S)-N-((2S,5S,8S,11S,14S,17S,20S,23S)-20-((1H-imidazol-4-yl)methyl)-14-((1H-indol-3-yl)methyl)-23-(2-amino-2-oxoethyl)-11-(4-aminobutyl)-17-((R)-1-hydroxyethyl)-5-(hydroxymethyl)-2,8-dimethyl-1,4,7,10,13,16,19,24,25,28-decaoxo-3,6,9-,12,15,18,21,22,27-nonaazanonacosan-29-yl)-4-(hydroxyamino)-7-((S)-2-(hydroxyamino)-3-phenylpropanamido)-9-(methylthio)-5,6-dioxononanamide, acetate

H—Ala—Ser—Ala—Lys—Trp—Thr—His—Asn—Gly—Gly—

Glu—Met—Phe—OH

• XCH₃CO₂H

Synonyms: FGF7 (152-164) (human), Fibroblast Growth Factor 7 (152-164) (human)

Peptide Sequence: ASAKWTHNGGEMF-OH

MF: C₆₃H₉₀N₁₈O₁₉S • XC₂H₄O₂

FW: 1,435.6

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

Oligopeptide-41 (acetate) is supplied as a solid. A stock solution may be made by dissolving the oligopeptide-41 (acetate) in the solvent of choice, which should be purged with an inert gas. Oligopeptide-41 (acetate) is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Oligopeptide-41 is a cosmeceutical peptide.¹ It corresponds to amino acids 152-164 of human FGF7.² Formulations containing oligopeptide-41 have been used in hair care products.

References

1. Oshimura, E. and Sakamoto, K. Amino acids, peptides, and proteins. *Cosmetic science and technology: Theoretical principles and applications* 285-303 (2017).
2. Rudd, T.R., Preston, M.D., and Yates, E.A. The nature of the conserved basic amino acid sequences found among 437 heparin binding proteins determined by network analysis. *Mol. Biosyst.* **13**(5), 852-865 (2017).

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, 06/10/2026

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