

PRODUCT INFORMATION



Klotho-derived Peptide 6 (acetate)

Item No. 41263

Formal Name: L-glutaminy-L-prolyl-L-valyl-L-valyl-L-threonyl-L-leucyl-L-tyrosyl-L-histidyl-L-tryptophyl-L- α -aspartyl-L-leucyl-L-prolyl-L-glutaminy-L-arginyl-L-leucyl-L-glutaminy-L- α -aspartyl-L-alanyl-L-tyrosylglycylglycyl-L-tryptophyl-L-alanyl-L-asparaginy-L-arginyl-L-alanyl-L-leucyl-L-alanyl-L- α -aspartyl-L-histidine, acetate
H—Gln—Pro—Val—Val—Thr—Leu—Tyr—His—Trp—Asp—Leu—Pro—Gln—Arg—Leu—Gln—Asp—Ala—Tyr—Gly—Gly—Trp—Ala—Asn—Arg—Ala—Leu—Ala—Asp—His—OH
• XCH₃CO₂H

Synonyms: Klotho (186-215) (human), KP6

Peptide Sequence: QPVVTLYHWDLPQRLQDAYGGWANRALADH-OH

MF: C₁₅₉H₂₃₂N₄₆O₄₄ • XC₂H₄O₂

FW: 3,491.8

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.

Description

Klotho-derived peptide 6 (KP6) is a synthetic peptide fragment that corresponds to amino acids 186-215 of the anti-aging protein Klotho.¹ It inhibits Wnt1-induced increases in fibronectin and α -smooth muscle actin (α -SMA) levels in HKC-8 human proximal tubular cells. KP6 (5 μ g/ml) also inhibits Wnt3a-induced activation of β -catenin in a reporter assay using HEK293T cells. *In vivo*, KP6 (1 mg/kg) increases survival and decreases urine albumin levels, renal hypertrophy, and glomerular injury in a mouse model of diabetic kidney disease induced by streptozotocin (STZ; Item No. 13104), advanced oxidation protein products, and unilateral nephrectomy.

Reference

1. Chen, X., Tan, H., Xu, J., *et al.* Klotho-derived peptide 6 ameliorates diabetic kidney disease by targeting Wnt/ β -catenin signaling. *Kidney Int.* **102**(3), 506-520 (2022).

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, 07/16/2025

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