

PRODUCT INFORMATION



BMP2-derived Peptide (trifluoroacetate salt)

Item No. 36948

Synonyms:	Bone Morphogenic Protein 2-derived Peptide, KIPKASSVPTELSAISTLYL-NH ₂	H-Lys-Ile-Pro-Lys-Ala-Ser-Ser-Val-Pro-Thr- Glu-Leu-Ser-Ala-Ile-Ser-Thr-Leu-Tyr-Leu- Gly-Gly-Cys-NH ₂
MF:	C ₁₀₄ H ₁₇₆ N ₂₆ O ₃₂ S • XCF ₃ COOH	
FW:	2,334.8	
Purity:	≥95%	
Supplied as:	A solid	
Storage:	-20°C	• XCF ₃ COOH
Stability:	≥4 years	

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

BMP2-derived peptide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the BMP2-derived peptide (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

BMP2-derived peptide is a synthetic peptide that corresponds to residues 73-92 of bone morphogenetic protein 2 (BMP2) and has a role in osteogenesis.¹ It increases alkaline phosphatase (ALP) activity in C3H/10T1/2 murine mesenchymal cells when used at concentrations of 100 and 200 µg/well. BMP2-derived peptide (100 µg/well) increases mRNA expression of the gene encoding osteocalcin in C3H/10T1/2 cells. It induces osteoblast-like cell differentiation and ectopic calcification when conjugated to an alginate gel in rats when administered via a submuscular implant.²

References

1. Saito, A., Suzuki, Y., Ogata, S.I., *et al.* Activation of osteo-progenitor cells by a novel synthetic peptide derived from the bone morphogenetic protein-2 knuckle epitope. *Biochim. Biophys. Acta* **1651**(1-2), 60-67 (2003).
2. Saito, A., Suzuki, Y., Ogata, S.-I., *et al.* Prolonged ectopic calcification induced by BMP-2-derived synthetic peptide. *J. Biomed. Mater. Res. A* **70**(1), 115-121 (2004).

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, 12/21/2022

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