

# BMP-4 (human), (recombinant)

Bone morphogenetic protein 4 (BMP-4) is a member of the bone morphogenetic growth factor family (BMP) and functions as a critical component in the culture and differentiation of stem cells. Active BMP-4 forms a homodimer or a heterodimer with related BMPs. BMP-4 is integral for the differentiation of mesoderm and ectoderm from pluripotent stem cells (iPSCs and ESCs) in stem cell cultures. BMP-4 is also used in culturing cardiomyocytes, early hematopoietic cells, hepatocytes, bone, cartilage, adipocytes, and lung cells, as well as inhibiting early neuronal differentiation.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-PRT259-0020	20µg
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)

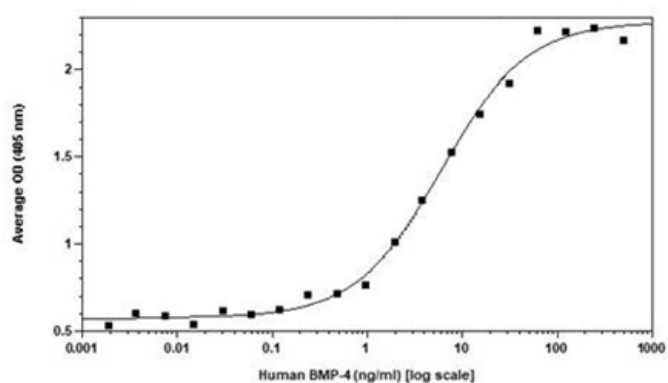


Figure 1. Human BMP-4 induced alkaline phosphatase activity in ATDC-5 cells.

## Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA solution.
Long Term Storage	-80°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Bone morphogenetic protein 4
Endotoxin Content	≤ 1 EU/μg
Formulation	Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing 0.1% Trifluoroacetic Acid (TFA).
MW	Dimer, 13.1/26.2 kDa (116/232 amino acids).
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in genetically modified CHO cells.
UniProt ID	P12644

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)