

Bone morphogenetic protein 4 (human), (recombinant)

Authentic glycosylation

BMP4 is a member of the bone morphogenetic protein family which is part of the TGF- β superfamily. The superfamily includes large families of growth and differentiation factors. Bone morphogenetic proteins were originally identified by an ability of demineralized bone extract to induce endochondral osteogenesis *in vivo* in an extraskeletal site. BMP4 plays an important role in the onset of endochondral bone formation in humans, and a reduction in expression has been associated with a variety of bone diseases, including the heritable disorder Fibrodysplasia Ossificans Progressiva.

Ordering Information

[Order Online »](#)

ENZ-PRT201-0010	10 μ g
-----------------	------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	The reconstituted solution can be stored at +4°C for 1 week.
Handling	After reconstitution, prepare aliquots and store at -20°C. For long term storage it is recommended to add a carrier protein (0.1% BSA or serum).
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	BMP-4
Appearance	White powder.
Formulation	Lyophilized from 0.7 mg/mL in 2X PBS containing 6% ethanol.
MW	34kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute in sterile 4 mM HCl containing 0.1% endotoxin free recombinant HSA.
Source	Produced in HEK cells. Glycosylated disulfide linked homodimer.
UniProt ID	P12644



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-
fr@enzolifesciences.com](mailto:info-fr@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)