

Bone morphogenetic protein 2 (human), (recombinant)

Authentic glycosylation

BMP2 belongs to the transforming growth factor-beta (TGFB) superfamily or structurally related signaling proteins. Bone morphogenic protein induces bone and cartilage formation. BMP2 is a candidate gene for the autosomal dominant disease of fibrodysplasia (myositis) ossificans progressiva.

Ordering Information

[Order Online »](#)

ENZ-PRT200-0010	10µg
-----------------	------

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-2
Appearance	White powder.
Formulation	Lyophilized from 1 mg/mL in 1X PBS.
MW	30-38kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Sterile filtered.
Reconstitution	Reconstitute with sterile 4 mM HCl containing 0.1% endotoxin-free recombinant HSA.
Source	Produced in HEK cells. Glycosylated disulfide-linked homodimer.
Specific Activity	The specific activity was ≤60ng/mL as determined by the dose dependent induction of alkaline phosphatase production in the ATDC-5 cell line.

UniProt ID

P12643



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Last modified: May 29, 2024