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 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-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

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com