

Recombinant Human BMP-4

Bone morphogenetic protein 4 (BMP-4) is a growth factor protein that is part of the transforming growth factor-beta superfamily. Responsible for the onset of endochondral bone formation, as well as muscle development, bone mineralization, mesoderm induction, tooth development, limb formation, fracture repair, and ureteric bud development. Dysregulation has been associated with a variety of bone diseases.

Materials Provided

Catalog Number:	92-1267
Quantity Supplied:	1 vial(s), 5 µg per vial
Reconstitution Buffer:	1 mL

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	31-36 kg/mol
Source:	Recombinant
Purity:	≥ 97%
Endotoxin Level:	< 1EU/µg
Formulation:	N/A
Storage:	-20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	100 µg/mL
Reconstitution Volume:	0.05 mL
Reconstitution Solvent:	Reconstitution Buffer

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com