

# PRODUCT INFORMATION



## BMP2 (recombinant)

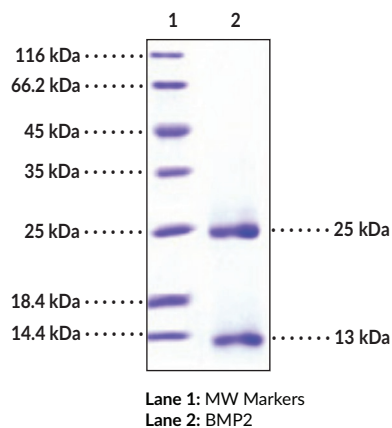
Item No. 32050

### Overview and Properties

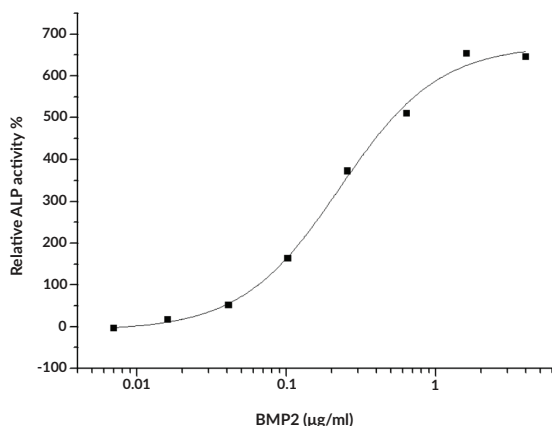
**Synonym:** Bone Morphogenic Protein 2  
**Source:** Active recombinant BMP2 expressed in *E. coli*  
**Amino Acids:** 283-396  
**Uniprot No.:** P12643  
**Molecular Weight:** 13 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥1 year  
**Purity:** ≥95% estimated by SDS-PAGE  
**Supplied in:** Lyophilized from sterile 30% CAN and 0.1% TFA, pH 2.9  
**Bioactivity:** See figures for details

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Images



**SDS-PAGE Analysis of BMP2.** This protein has a calculated molecular weight of 13 kDa. By SDS-PAGE, under non-reducing conditions, the protein has apparent molecular weights of approximately 13 and 25 kDa corresponding to the monomer and dimer, respectively.



**ALP Induction by BMP2.** BMP2 is measured by its ability to induce alkaline phosphatase production by MC3T3-E1 mouse osteoblastic cells. The  $EC_{50}$  value for this effect is typically 0.1-0.5 µg/ml.

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/15/2020

**CAYMAN CHEMICAL**

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

Bone morphogenetic protein 2 (BMP2) is a member of the TGF- $\beta$  superfamily of growth factors that has roles in osteogenesis and embryogenesis.<sup>1</sup> It is composed of an N-terminal prodomain that regulates BMP2 folding and secretion and a C-terminal mature domain that contains the receptor binding region. BMP2 is expressed by chondrocytes, osteocytes, osteoblasts, and vascular endothelial cells and is synthesized in the cytosol as a proprotein dimer.<sup>2,3</sup> Cleavage of the proprotein dimer by proprotein convertase subtilisin kexins (PCSKs) generates the N- and C-terminal domains, which remain non-covalently associated and homodimerize or heterodimerize with BMP4 or BMP7 prior to secretion.<sup>2</sup> BMP2 preferentially binds to BMP type 1 receptors, also known as activin receptor-like kinases (ALKs), which are expressed by mesenchymal or hematopoietic stem cells, chondrocytes, osteoblasts, osteoclasts, and myoblasts, and activates SMAD-dependent and -independent gene transcription.<sup>1,4</sup> BMP2 induces production of collagen type I  $\alpha$  and alkaline phosphatase, a marker of bone formation, as well as induces differentiation and proliferation of bone marrow-derived mesenchymal stem cells *in vitro*.<sup>5</sup> Genome-wide deletion of *Bmp2* in mice induces amnion, chorion, and cardiac developmental defects and is embryonic lethal.<sup>1</sup> Limb-specific *Bmp2* conditional knockout mice exhibit spontaneous fractures and impaired fracture healing.<sup>6</sup> Serum and synovial fluid BMP2 levels are increased in patients with osteoarthritis and are positively correlated with disease severity.<sup>7</sup> Formulations containing BMP2 have been used in treatment of long bone fractures. Cayman's BMP2 (recombinant) protein can be used for cell-based assay applications. This protein consists of 115 amino acids that correspond to the mature form of human, canine, mouse, rat, and rhesus BMP2 and has a calculated molecular weight of 13 kDa. By SDS-PAGE, under non-reducing conditions, the protein has apparent molecular weights of approximately 13 and 25 kDa, corresponding to the monomer and dimer, respectively.

## References

---

1. Wang, R.N., Green, J., Wang, Z., *et al.* Bone morphogenetic protein (BMP) signaling in development and human diseases. *Genes Dis.* **1**(1), 87-105 (2014).
2. McBride-Gagy, S., McKenzie, J.A., Buettmann, E.G., *et al.* Bmp2 conditional knockout in osteoblasts and endothelial cells does not impair bone formation after injury or mechanical loading in adult mice. *Bone* **81**, 533-543 (2015).
3. Constam, D.B. Regulation of TGF $\beta$  and related signals by precursor processing. *Semin. Cell Dev. Biol.* **32**, 85-97 (2014).
4. Lin, S., Svoboda, K.K.H., Feng, J.Q., *et al.* The biological function of type I receptors of bone morphogenetic protein in bone. *Bone Res.* **4**, 16005 (2016).
5. Lysdahl, H., Baatrup, A., Foldager, C.B., *et al.* Preconditioning human mesenchymal stem cells with a low concentration of BMP2 stimulates proliferation and osteogenic differentiation *in vitro*. *Biores. Open Access* **3**(6), 278-285 (2014).
6. Tsuji, K., Bandyopadhyay, A., Harfe, B.D., *et al.* BMP2 activity, although dispensable for bone formation, is required for the initiation of fracture healing. *Nat. Genet.* **38**(12), 1424-1429 (2006).
7. Liu, Y., Hou, R., Yin, R., *et al.* Correlation of bone morphogenetic protein-2 levels in serum and synovial fluid with disease severity of knee osteoarthritis. *Med. Sci. Monit.* **21**, 363-370 (2015).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM