

# Mouse FGF basic (recombinant)



**Alias:** FGF-2 **Catalog #:** 6506

**Size:** 5 ug **Research Use Only**

**Molecular Weight:** 16.2 kDa

**Source:** Yeast. Recombinant Mouse FGF basic (FGF-2) was produced in yeast and, therefore, does not have endotoxin. It is naturally folded and post-translationally modified.

**Formulation:** Lyophilized without carrier protein.

**Purity:** >95% as visualized by SDS-PAGE analysis.

**Purification:** Ion-exchange chromatography.

**Bioactivity:** In testing

**Entrez Gene ID:** 14173

**Number of Amino Acids:** 144

**Amino Acid Sequence:** ALPEDGGAAP PPGHFKDPKR LYCKNGGFFL RIHPDGRVDG VREKSDPHVK LQLQAEERGV  
VSIKGVCANR YLAMKEDGRL LASKCVTEEC FFERLESNN YNTYRSRKYS SWYVALKRTG  
QYKLGSKTGP GQKAILFLPM SAKS (144)

**Country of Origin:** USA

**Reconstitution:** Reconstitute with sterile phosphate-buffered saline containing at least 0.1% carrier protein.

**Stability and Storage:** Stable for up to twelve months from date of receipt at -20°C. Stable for at least 3 months when stored in working aliquots with a carrier protein at -20°C. Avoid repeated freeze/thaw cycles.

**Applications:** The FGF basic protein can be used in cell culture, as a FGF basic ELISA Standard, and as a Western Blot Control.

**Background:** Fibroblast growth factors, or FGFs, are a family of growth factors involved in angiogenesis, wound healing, and embryonic development. The FGF family members are heparin-binding proteins and interactions with cell-surface-associated heparan sulfate proteoglycans have been shown to be essential for FGF signal transduction. There are currently 24 members of the FGF family. FGF basic (FGF2) and FGF acidic (FGF1) are multipotential factors that stimulate and support proliferation, migration and differentiation.

F17-6358-4-A Data Sheet; Effective: 1/10/14; Supersedes: None; Page 1 of 1; Recombinant Mouse FGF basic (FGF-2) updated on: 1/29/2014