

Mouse IFN beta (recombinant)



Alias: none **Catalog #:** 6541

Size: 5 ug **Research Use Only**

Molecular Weight: 19.7 kDa

Source: Yeast. Recombinant Mouse IFN beta 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: 15977

Number of Amino Acids: 161

Amino Acid Sequence: INYKQLQLQE RTNIRKCQEL LEQLNGKINL TYRADFKIPM EMTEKMQKSY TAFAIQEMLQ
NVFLVFRNNF SSTGWNETIV VRLDELHQQ TVFLKTVLEE KQEERLTWEM SSTALHLKSY
YWRVQRYLKL MKYNSYAWMV VRAEIFRNFL IIRRLTRNFQ N (161)

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 IFN beta protein can be used in cell culture, as a IFN beta ELISA Standard, and as a Western Blot Control.

Background: Type I interferons (IFN alpha, IFN beta) belong to the helical cytokine superfamily, which includes growth hormones, interleukins, several colony-stimulating factors plus other regulatory molecules. All function as regulators of cellular activity by interacting with cell-surface receptors and activating various signaling pathways. IFN beta produces antiviral, antibacterial, and anticancer properties. RANKL (TNFSF11) has been shown to induce the production of IFN beta.

F17-6358-4-A Data Sheet; Effective: 1/10/14; Supersedes: None; Page 1 of 1; Recombinant Mouse IFN beta updated on: 1/29/2014