

Bovine BAFF (recombinant)



Alias: TNFSF13B **Catalog #:** 6536

Size: 5 ug **Research Use Only**

Molecular Weight: 17.1 kDa

Source: Yeast. Recombinant Bovine BAFF (TNFSF13B) 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: 504507

Number of Amino Acids: 152

Amino Acid Sequence: ALQEAETVT QDCLQLIADS DTPTIRKGAY TFVPWLLSFK RGRALEEKEN KILVKETGYF
FIYGQVLYTD NTFAMGHLIQ RKKVHVFGDE LSLVTLFRCI QNMPETLPNN SCYSAGIAKL
EEGDELQLAI PREDAKISRD GDGTFFGALK LL (152)

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 Bovine BAFF protein can be used in cell culture, as an BAFF ELISA Standard, and as a Western Blot Control.

Background: BAFF (B-cell Activation Factor), also known as Tumor Necrosis Factor Superfamily member 13B (TNFSF13B), is a member of the tumor necrosis factor (TNF) ligand family. Nineteen cytokines have been identified as part of the TNF family on the basis of sequence, functional, and structural similarities. Family members include TNF beta (TNFSF1), TNF alpha (TNFSF2), Lymphotoxin beta (TNFSF3), OX40 Ligand (TNFSF4), CD40 Ligand (TNFSF5), Fas Ligand (TNFSF6), CD27 Ligand (TNFSF7), CD30 Ligand (TNFSF8), 4-1BB Ligand (TNFSF9), TRAIL (TNFSF10), TRANCE/RANKL (TNFSF11), TWEAK (TNFSF12), APRIL (TNFSF13), BAFF (TNFSF13B), LIGHT (TNFSF14), TL1A/VEGI (TNFSF15), and GITR Ligand (TNFSF18). BAFF is expressed in B cell lineage cells and acts as a potent B cell activator. It has been also shown to play an important role in the proliferation and differentiation of B cells.

F17-6358-4-A Data Sheet; Effective: 1/10/14; Supersedes: None; Page 1 of 1; Recombinant Bovine BAFF (TNFSF13B) updated on: 1/29/2014