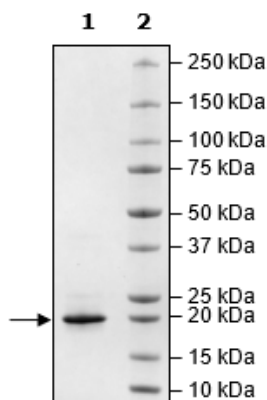


## Product Information

<b>Description:</b>	Recombinant human BAFF (B-cell-activating factor, also known as TNFSF13B, Tumor necrosis factor ligand superfamily member 13B), encompassing amino acids 134-285. This construct contains an N-terminal His-tag (6xHis) followed by an Avi-Tag™. The recombinant protein was affinity purified.
<b>Background:</b>	B-cell activating factor (BAFF) is a cytokine regulating B cell development, survival, and antibody production. It is a key player in the immune response and has implications in autoimmune diseases. It is a ligand for receptors TNFRSF13B/TACI, TNFRSF17/BCMA, and TNFRSF13C/BAFFR. BAFF normally maintains immune balance, and overproduction of BAFF leads to various auto-immune diseases. It is, therefore, a therapeutic target. Drugs that inhibit BAFF, such as belimumab, have been approved for the treatment of certain autoimmune disorders.
<b>Species:</b>	Human
<b>Construct:</b>	BAFF (His-Avi-134-285(end))
<b>Concentration:</b>	0.82 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>MW:</b>	20 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Genbank Accession:</b>	NM_006573
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Assay Conditions:</b>	Assay was performed according to BAFF:BCMA[Biotinylated] Inhibitor Screening Assay Kit (BPS Bioscience #79667) with BCMA titrated at various concentrations.
<b>Applications:</b>	Useful for binding studies and screening inhibitors of ligand binding to its receptor.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



BAFF:BCMA[Biotinylated] Binding

