

Recombinant TRβ1 protein

Catalog No: 31139

Expressed In: *E. coli*

Quantity: 10 µg

Concentration: 0.15 µg/µl

Source: Human

Buffer Contents: 10 µg of Recombinant TRβ1 protein in Dilution Buffer AM1 (20 mM Tris-Cl (pH 8), 20% glycerol, 100 mM NaCl, 1 mM DTT and 0.2 mM EDTA).

Protein Details: Recombinant TRβ1 protein is isolated from an *E. coli* strain that carries the coding sequence of the human TRβ1 isoform under the control of a T7 promoter (accession number NM 000461) and purified by a combination of affinity and gel filtration chromatography. TRβ1 protein has an N-terminal His-Tag and is purified free of contaminating proteases, endonucleases and exonucleases, and is visually determined to be > 95% pure on an SDS-polyacrylamide gel.

Application Notes: Recombinant TRβ1 is suitable for *in vitro* transcription, chromatin remodeling and protein-protein interaction assays. 20 ng is sufficient for reconstituted transcription assays and 100 ng is sufficient for protein-protein interaction studies. The molecular weight of the protein is ~45 kDa.

NOTE: The presence of Poly [d(I-C)] in buffers may affect protein functionality and should be avoided.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.