

## Recombinant APEX protein

---

**Catalog No:** 31245

**Expressed In:** *E. coli*

**Quantity:** 10 µg

**Concentration:** 0.2 µg/µl

**Source:** Human

**Buffer Contents:** 10 µg of recombinant protein supplied at a concentration of 0.2 µg/µl in 20 mM Tris-HCl, pH 7.3, 300 mM KCl, 0.2 mM EDTA and 20% glycerol.

**Background:** APEX, APE/Ref-1 apurinic/aprimidinic endonuclease (also known as APE, Ref1, AP endonuclease I or APEX nuclease) is a DNA repair enzyme that removes the blocking group from the 3'-termini of the DNA strand breaks generated by ionizing radiation. The apurinic/aprimidinic (AP) sites can prevent normal DNA replication.

**Protein Details:** Recombinant APEX protein was expressed in *E. coli* as the full-length human protein (accession number AAB22977) with an N-terminal FLAG tag. The molecular weight of the protein is ~38 kDa.

**Application Notes:** Recombinant APEX protein is suitable for use in DNA repair assays. We recommended to start with a range of protein concentrations (*e.g.* 5 ng, 50 ng, and 500 ng) to determine the optimal amount needed for each assay.

**Storage and Guarantee:** This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.