

Recombinant DNA Topoisomerase I protein

Catalog No: 31239

Expressed In: Baculovirus

Quantity: 10 µg

Concentration: 0.15 µg/µl

Source: *Drosophila melanogaster*

Buffer Contents: 10 µg of recombinant protein supplied at a concentration of 0.15 µg/µl in 20 mM Tris-HCl, pH 7.3, 125 mM NaCl, 0.2 mM EDTA and 20% glycerol.

Background: DNA topoisomerase I (TOP1) is an enzyme involved in the state of DNA during transcription. It catalyzes the breaking and rejoining of single stranded DNA, thus altering the topology of DNA.

Protein Details: Recombinant DNA Topoisomerase I protein was expressed in Sf9 insect cells as the full-length *Drosophila melanogaster* protein (accession number AAA28951) with an N-terminal FLAG tag. The molecular weight of the protein is ~135 kDa.

Application Notes: Recombinant DNA Topoisomerase I protein is suitable for use in assays that study DNA and chromatin topological changes. A recommended starting point is 5-20 ng recombinant protein per reaction.

Storage and Guarantee: 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.