

Recombinant YY1 protein

Catalog No: 31332

Source: Human

Quantity: 10 µg

Expressed In: *E. coli*

Buffer: The protein is supplied at a concentration of 0.25 µg/µl in 20 mM Tris-HCl (pH7.3), 300 mM KCl, 0.2 mM EDTA and 20% glycerol.

Contents: 10 µg recombinant YY1 protein.

Protein Details:

YY1 (Yin Yang 1) is a transcription factor that belongs to the GLI-Kruppel class of zinc finger transcription factors and is ubiquitously expressed. YY1 is known to have a fundamental role in normal biologic processes such as embryogenesis, differentiation, replication and cellular proliferation. YY1 exerts its effects on genes involved in these processes via its ability to initiate, activate or repress transcription depending upon the context in which it binds. Since expression and function of YY1 are known to be intimately associated with progression through phases of the cell cycle, the physiologic significance of YY1 activity has been applied to models of tumor biology. It can repress and activate a number of gene promoters and direct histone deacetylases (HDACs) and histone acetyltransferases (HATs). YY1 is known to facilitate silencing by the Polycomb protein complex.

Dilution of Protein:

Recombinant YY1 is supplied in solution and should be kept as a solution in order to maintain full activity.

Application:

Recombinant YY1 protein is suitable for use in *in vitro* transcription assays. We recommend 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 *in vitro* transcription assay.

Storage:

Recombinant proteins in solution are temperature sensitive and should be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage.

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.