

Recombinant SOX9 protein



Catalog No: 82079, 82679

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Source: Human

Buffer Contents: Recombinant SOX9 protein is supplied in 25 mM HEPES pH 7.5, 300 mM NaCl, 20% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

Background: **SOX9** is a transcription factor that plays a key role in chondrocytes differentiation and skeletal development. Specifically binds the 5'-ACAAAG-3' DNA motif present in enhancers and super-enhancers and promotes expression of genes important for chondrogenesis, including cartilage matrix protein-coding genes COL2A1, COL4A2, COL9A1, COL11A2 and ACAN, SOX5 and SOX6.

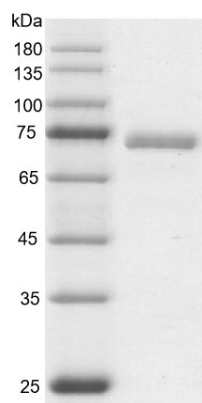
Protein Details: Recombinant SOX9 protein was expressed in a baculovirus expression system as the full length protein (accession number NP_000337.1) with an N-terminal FLAG tag. The molecular weight of SOX9 is 57.41 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for this product is shown.

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.

SOX9



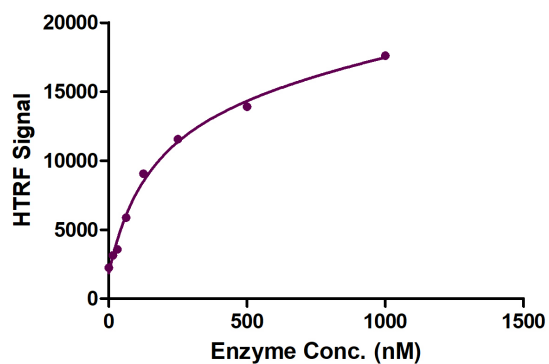
Recombinant SOX9 protein

10% SDS-PAGE Coomassie staining

MW: 57.41 kDa,

Purity: $\geq 95\%$

SOX9 Titration



HTRF Assay for Recombinant SOX9 activity

1 μM substrate DNA (5' biotin-CGACCATTGTTTCAG 3') was incubated with different concentrations of SOX9 protein in reaction buffer containing 50 mM HEPES NaOH pH 7.0, 0.1% BSA at room temperature for 1 hour. Anti-FLAG antibody was used to detect reaction products.