



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-44460

NBP1-44460 Protocol

SDS-Page protocol for Sorbitol Dehydrogenase Recombinant Protein (NBP1-44460)

SDS-PAGE (NBP1-44460): https://www.novusbio.com/products/sorbitol-dehydrogenase-recombinant-protein_nbp1-44460

1. Prepare a 1.5 ml assay buffer. The final concentrations are 78 mM Triethanolamine, 183 mM D-Fructose, 0.2 mM beta-nicotinamide dinucleotide and 0.003 % (w/v) bovine serum albumin. <p/>
2. Add 50 ul of recombinant SORD protein in various concentrations (1 ug, 2 ug, 5 ug) in assay buffer. <p/>
3. Mix by inversion and record the decrease at A340nm for 5 minutes.