



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-40395

NBP1-40395 Protocol

SDS-Page protocol for Aldolase A Biologically Active Protein (NBP1-40395)

SDS-PAGE (NBP1-40395): https://www.novusbio.com/products/aldolase-a-biologically-active-protein_nbp1-40395

1. Prepare a 1.45 ml reaction mixture with the following concentrations: 45 mM Tris pH8.0, 0.95mM fructose 1,6-diphosphate, 0.065 mM beta-nicotinamide adenine dinucleotide, 5 unit alpha-glycerophosphate dehydrogenase/triosephosphate isomerase.

2. Add 50 ul of recombinant Aldolase A protein with 1ug, 2ug in reaction mixture.

3. Record the decrease in A340nm for 5 minutes at 25C.