



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-30262

NBP1-30262 Protocol

SDS-Page protocol for Enolase 1 Recombinant Protein (NBP1-30262)

SDS-PAGE (NBP1-30262): https://www.novusbio.com/products/enolase-1-recombinant-protein_nbp1-30262

1. Prepare 1,450ul assay buffer. The final concentrations are 81 mM Triethanolamine, 1.9 mM 2-phosphoglycerate, 0.12 mM beta- NADPH, 25 mM magnesium sulfate, 100 mM potassium chloride, 1.3 mM ADP, 4 unit pyruvate kinase, 6 unit L-lactic dehydrogenas. <p/>

2. Add 50ul of recombinant protein alpha-enolase in various concentrations(0.25 ug, 0.5 ug, and 0.1 ug) in assay buffer. <p/>

3. Mix by inversion and load 200 ul of reaction mix in to a plate well. <p/>

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