



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-99038

NBP1-99038 Protocol

SDS-Page protocol for MDH1 Recombinant Protein (NBP1-99038)

SDS-PAGE (NBP1-99038): https://www.novusbio.com/products/mdh1-recombinant-protein_nbp1-99038

1. Prepare a 1.45 ml assay buffer. The final concentrations are 100 mM Potassium phosphate, 0.13mM beta-nicotinamide adenine dinucleotide, reduced form, 0.25mM oxalacetic acid. <p/>
2. Add 50 ul of recombinant MDH protein with various concentrations (0.015 ug, 0.03 ug) in assay buffer. <p/>
3. Mix by inversion and record the decrease at A340nm for approximately 5 minutes.