



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-72281

NBP1-72281 Protocol

SDS-Page protocol for Dihydrofolate Reductase/DHFR Biologically Active Protein (NBP1-72281)

SDS-PAGE (NBP1-72281): https://www.novusbio.com/products/dihydrofolate-reductase-dhfr-biologically-active-protein_nbp1-72281

1. Prepare a 1.55 ml assay buffer. The final concentrations are 50 mM potassium phosphate, 0.072 mM dihydrofolic acid, 0.1 mM beta-nicotinamide dinucleotide phosphate and 0.003 % (w/v) bovine serum albumin. <p/>
2. Add 50 ul of recombinant DHFR protein in various concentrations (2 ug, 5 ug) in assay buffer. <p/>
3. Mix by inversion and record the decrease at A340nm for 5 minutes.