



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-78817

NBP1-78817 Protocol

SDS-PAGE (NBP1-78817)

1. Prepare a 1,450 ul assay buffer. The final concentrations are 75mM Potassium phosphate (pH 7.5), 2.6mM EDTA 1mM glutathione, 0.09mM beta-NADPH, 0.13% BSA.

2. Add 50 ul of recombinant GSR protein with 0.018 ug, 0.037 ug and 0.075 ug in assay buffer

3. Mix and load 200 ul of reaction mix in to a plate well.

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