



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-50895

NBP1-50895 Protocol

SDS-Page protocol for PYCR1 Protein (NBP1-50895)

SDS-PAGE (NBP1-50895): https://www.novusbio.com/products/pycr1-protein_nbp1-50895

1. Prepare an assay buffer with the following concentrations:

200 mM Tris-HCl (pH10.0) and 10 mM NAD, 20mM L-proline, and load 150 ul into each well. <p/>

2. Add 50 ul of 1.25 ug and 2.5 ug recombinant PYCR1 protein to each well. <p/>

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