



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-30292

NBP1-30292 Protocol

SDS-Page protocol for PKM2 Recombinant Protein (NBP1-30292)

SDS-PAGE (NBP1-30292): https://www.novusbio.com/products/pkm2-recombinant-protein_nbp1-30292

1. Prepare a 150ul reaction mixture into a suitable container.

- Reaction mixture: 72mM Imidazole (pH 7.5), 7.6mM ADP, 15mM MgCl₂, 72mM KCl, 0.2mM Beta-NADH.

2. Incubate for 10 min at 37C.

3. Add 50ul of 20 units/ml LDH, 50ul of 5mM PEP, and incubate for 10 min at 37C.

4. Add 50ul of recombinant PKM2 protein (in 0.1% BSA / PBS) with various concentrations (0.03ug -1ug).

5. Read the decrease in A340nm in kinetic mode for 10 minutes. <p/>