



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP1-49306

NBP1-49306 Protocol

SDS-PAGE (NBP1-49306)

1. Prepare 150 μ l assay buffer. The final concentrations are 50 mM Tris-HCl (pH 7.5), 100 μ M cytidine.

 2. Add 50 μ l recombinant CDA protein with concentration 0.25 μ g in assay buffer.

 3. Mix by inversion and records the decrease at A280nm for approximately 5 minutes. <p/>
- Cytidine millimolar extinction coefficient: 5.27 (mM-1.cm-1) 100 μ M / 50 mM Tris-HCl (pH 7.5)