



Orders: [orders@novusbio.com](mailto:orders@novusbio.com)

Support: [technical@novusbio.com](mailto:technical@novusbio.com)

Web: [www.novusbio.com](http://www.novusbio.com)

Protocols, Publications, Related Products, Reviews and more:

[www.novusbio.com/NBP1-98907](http://www.novusbio.com/NBP1-98907)

## NBP1-98907 Protocol

### SDS-Page protocol for Catalase Recombinant Protein (NBP1-98907)

---

SDS-PAGE (NBP1-98907): [https://www.novusbio.com/products/catalase-recombinant-protein\\_nbp1-98907](https://www.novusbio.com/products/catalase-recombinant-protein_nbp1-98907)

1. Prepare an assay buffer with the following concentrations: 50 mM Potassium Phosphate (pH 8.0) and 0.15 % H<sub>2</sub>O<sub>2</sub>. <p/>
2. Load 150ul assay buffer into each well <p/>
3. Add 50ul of recombinant CAT protein with 0.25ug, 0.125ug to each well. <p/>
4. Record the decrease in A<sub>260nm</sub> for 5 minutes at 25C using UV plate.