



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/NBP1-72478](http://www.novusbio.com/NBP1-72478)

## NBP1-72478 Protocol

### SDS-PAGE (NBP1-72478)

1. Prepare a 200ul reaction mix into a suitable container: The final concentrations are 10mM Tris (pH7.5), 50mM NaCl, 2mM DTT, 1mM MnCl<sub>2</sub>, 10mM pNPP. <p/>
2. Pre-warm reaction mix in heat block at 37C.<p/>
3. Add recombinant DUSP3 protein solution with various concentrations (1ug, 2ug, 3ug) in 200ul reaction buffer. <p/>
4. Mix by inversion and incubate at 37C for 10minutes. <p/>
5. Stop the reaction by adding the 800ul of 0.1N NaOH and record the absorbance at 405nm.