



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/NBP1-45276](http://www.novusbio.com/NBP1-45276)

## NBP1-45276 Protocol

### SDS-Page protocol for UDP glucose dehydrogenase Biologically Active Protein (NBP1-45276)

---

SDS-PAGE (NBP1-45276): [https://www.novusbio.com/products/udp-glucose-dehydrogenase-biologically-active-protein\\_nbp1-45276](https://www.novusbio.com/products/udp-glucose-dehydrogenase-biologically-active-protein_nbp1-45276)

1. Prepare a 1470 ul assay buffer. The final concentrations are 93mM Tris-Hcl (pH 8.0), 0.2mM UDPG, and 1mM beta-NAD. <br>
2. Add 20 ul of recombinant protein UGDH with 1 ug, 2 ug and 5 ug in assay buffer<br>
3. Mix by inversion and load 200 ul of reaction mix in to a plate well.<br>
4. Record the decrease in A340nm for 5 minutes at 25C<br>