

## Data Sheet

### Peroxiredoxin 2 (PRDX2)

human, recombinant, N-terminal His-tag  
**Catalog #:** 70011  
**Lot#** 2003                      **Conc.** 2.1 mg/ml

**Formulated in:** 25 mM Tris-HCl, pH 8.0,  
 100 mM NaCl, 1.35 mM KCl, 0.05% Tween-  
 20, 50% glycerol, and 3 mM DTT.

**Stability:** >6 months at -80°C

**References:**

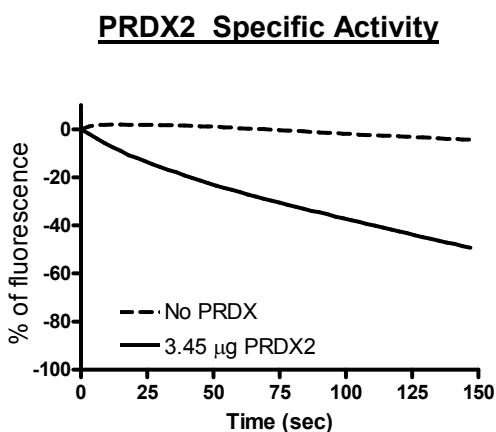
1. Peskin, A.V., *et al.*, *J. Biol. Chem.* **282** (16), 11885-11892 (2007)
2. Low, F.M., *et al.*, *Blood* **109** (6), 2611-2617 (2007)

**Description:** Human Peroxiredoxin 2, also known as PRDX2, full length (Gene bank accession # NM\_005809) with N-terminal His-tag, MW=22.7 kDa, expressed in Baculovirus infected Sf9 cell expression system.

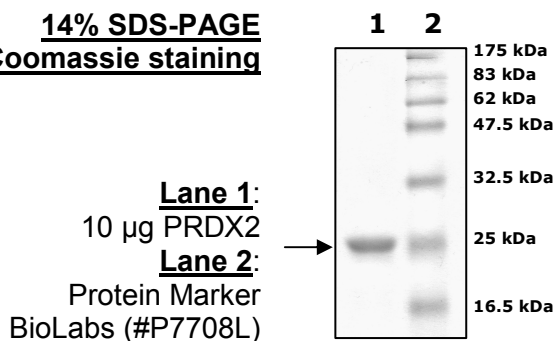
**Specific Activity:** 149 pmol/min/ug.  
 3.45 µg of PRDX2 was incubated in 50 mM HEPES (pH 7.0) containing 200 µM NADPH, 3 µM thioredoxin, and 1.5 µM thioredoxin reductase. The reaction mixture was incubated at 30°C for 5 min, followed by the addition of 0.22 mM H<sub>2</sub>O<sub>2</sub>. NADPH oxidation was monitored (fluorescence decrease) for the next 10 min.

**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance



**14% SDS-PAGE  
 Coomassie staining**



**MW:** 22.7 kDa  
**Purity:** >95%

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)