

Data Sheet

Peroxiredoxin I, His-tag

Human, recombinant, N-terminal His tag
Catalog #: 70010
Lot #: 160622 **Conc.:** 1.68 mg/ml

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

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Maksimov, V. et al. Genetika 42 (2), 274-277 (2006)
2. Jang, H.H et al. FEBS Lett. 580 (1), 351-355 (2006)

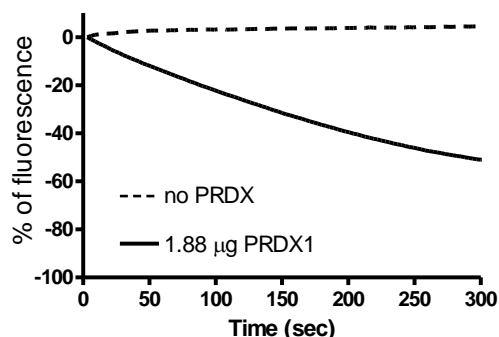
Description: Human Peroxiredoxin I (GenBank Accession No. NM_002574/NP_002565), full length, with N-terminal HIS6 tag, MW=23 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: >800 pmol/min/ug.
 1.88 μg of PRDX1 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_2O_2 . NADPH oxidation was monitored (fluorescence decrease) for the next 10 min.

Application: Useful for the study of biochemical function.

Quality Assurance

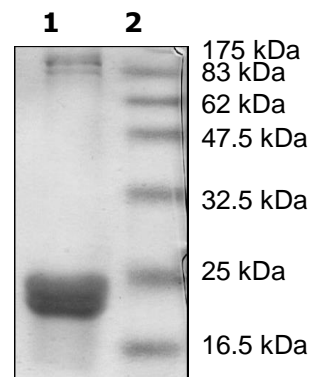
PRDX1 Specific Activity



**14% SDS-PAGE
Coomassie staining**

Lane 1:
20 μg Peroxiredoxin I
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 23 kDa.
Purity: $\geq 90\%$



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
 Please visit our website at: www.bpsbioscience.com