

Data Sheet

Peroxiredoxin I (K197R)

Human, recombinant, N-terminal His tag
Catalog #: 70021
Lot #: 111219 **Conc.:** 1.24 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 200 mM
Imidazole, 20% glycerol and 3 mM DTT.

Stability: >6 months at -80°C

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 (K197R), or PRDX1, GenBank Accession No. NM_002574/ NP_002565, a.a. 1-199 (end) full length protein containing a Lysine to Arginine mutation at position 197 and N-terminal His6 tag, MW=23 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Application:

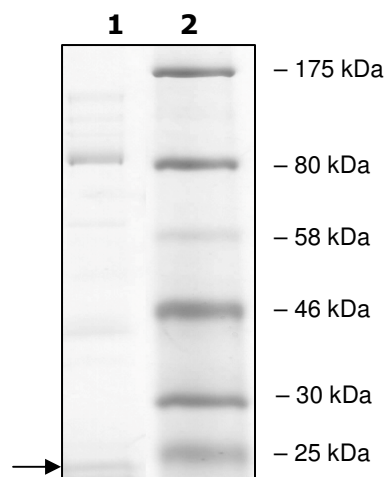
Useful for the study of biochemical function.

Quality Assurance

10% SDS-PAGE
Coomassie staining

Lane 1:
8.7 µg PRDX1 K197R
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 23 kDa
Purity: ≥92%



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