

NQO2

(QR2)

CATALOG NO.: NQO-11-312

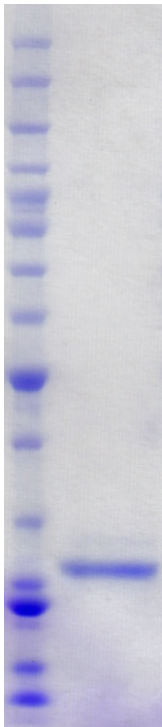
LOT NO.:

DESCRIPTION: Human recombinant NQO2 (NAD(P)H dehydrogenase quinone 2) protein (residues 1-231; Genbank Accession #NM_000904.4; MW = 26.0 kDa) expressed in *E. coli*.

PURITY: >90% by SDS-PAGE

SUPPLIED AS: $\mu\text{g}/\mu\text{L}$ in 50 mM Tris HCl, pH 7.5, 125 mM NaCl, 1 mM TCEP, 10 % glycerol

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry/ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 μl) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (12% acrylamide) of 4 μg of RBC NQO2. MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, **50**, 40, 30, 25, **20**, 15, 10 kDa.

This product is not intended for therapeutic or diagnostic use in animals or in humans.

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com