

## NQO1

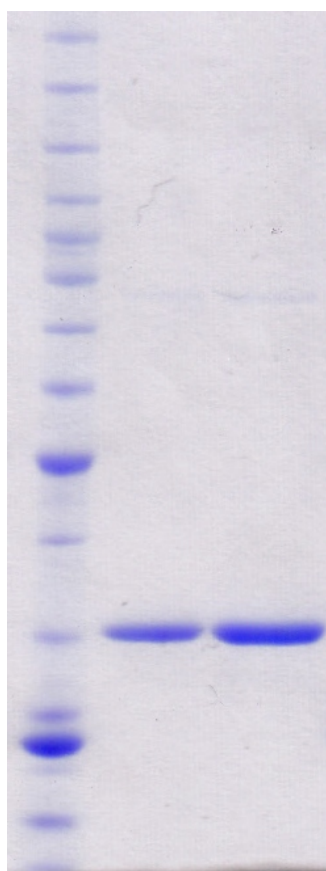
(QR1)

**CATALOG NO.:** NQO-11-314

**LOT NO.:**
**DESCRIPTION:** Human recombinant NQO1 (NAD(P)H dehydrogenase quinone 1) protein (residues 1-274; Genbank Accession # NM\_000903.2; MW = 31.0 kDa) expressed in *E. coli*.

**PURITY:** >90% by SDS-PAGE

**SUPPLIED AS:** \_ µg/µL in 50mM Tris-HCl, pH 7.5, 125mM NaCl, 10% glycerol, 1mM TCEP

**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 2 and 4 µg of RBC NQO1** 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