

NQO1 (QR1)

CATALOG NO.: NQO-11-314

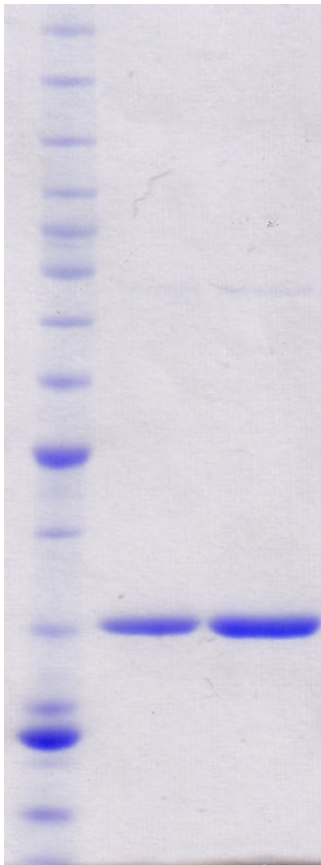
LOT NO.:

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: $_ \mu\text{g}/\mu\text{L}$ 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.