

NQO1 (His) (QR1)

CATALOG NO.: NQO-11-313

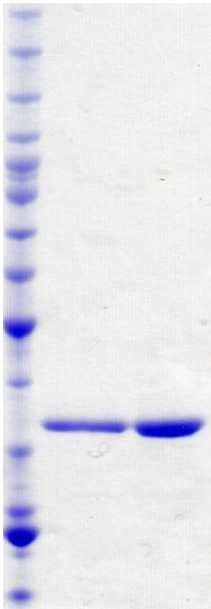
LOT NO.:

DESCRIPTION: Human recombinant NQO1 (NAD(P)H dehydrogenase quinone 1) protein (residues 1-274; Genbank Accession # NM_000903.2; MW = 33.5 kDa) expressed with an N-terminal His tag in *E. coli*.

PURITY: >95% 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 (His)
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.