

NAD(P)H dehydrogenase, quinone 2 (NQO2), transcript variant 3, (His)

**CATALOG NO.:** NQO-11-339

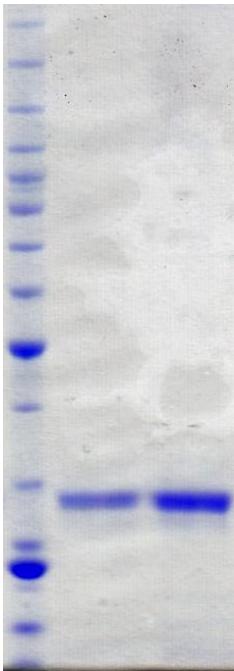
**LOT NO.:**

**DESCRIPTION:** Human recombinant NQO2 (residues 2-231; Genbank Accession # NM\_000904.4; MW = 28.5 kDa) expressed with an N-terminal His-tag in *E. coli*.

**PURITY:** >95% 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  $\mu\text{l}$ ) or storage of diluted enzyme is not recommended.



**Coomassie blue-stained SDS-PAGE (12% acrylamide) of 2 $\mu\text{g}$  and 4 $\mu\text{g}$  of RBC NQO2 (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.