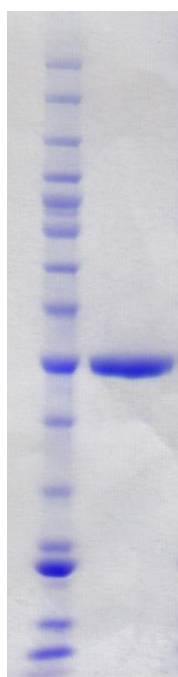


NQO2 (GST) (NAD(P)H dehydrogenase, quinone 2 (NQO2), transcript variant 3, (GST))

CATALOG NO.: NQO-11-340

LOT NO.:

DESCRIPTION: Human recombinant NQO2 (residues 2-231; Genbank Accession # NM_000904.4; MW = 52.8 kDa) expressed as an N-terminal GST-fusion protein 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 μl) or storage of diluted enzyme is not recommended.

Coomassie blue-stained SDS-PAGE (12% acrylamide) of 4 μg of RBC NQO2 (GST). 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