

Deoxyuridine Triphosphatase (DUT), transcript variant 1 (GST)

CATALOG NO.: NUC-11-335

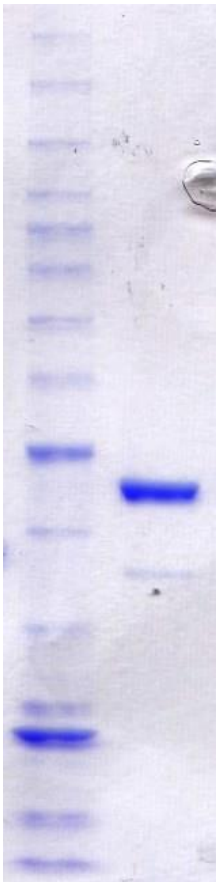
LOT NO.:

DESCRIPTION: Human recombinant DUT (residues 70-252; Genbank Accession #NM_001025248.1; MW = 46.2 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

PURITY: 90% by SDS-PAGE

SUPPLIED AS: $_ \mu\text{g}/\mu\text{L}$ in 50 mM Tris HCl, pH 7.5, 500 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 1 μg of RBC DUT (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.