

Nudix (Nucleoside Diphosphate linked moiety X)-type Motif 14 (NUDT14) (GST)

CATALOG NO.: NUC-11-343

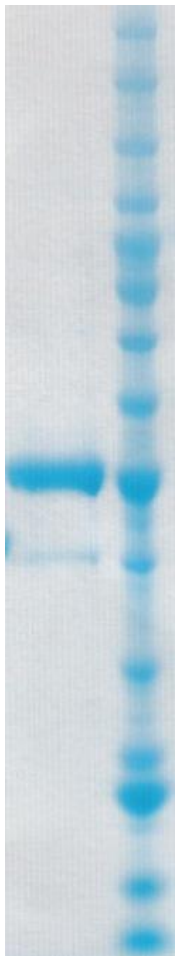
LOT NO.:

DESCRIPTION: Human recombinant NUDT14 (residues 2-222; Genbank Accession #NM_177533.4; MW = 50.9 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 2 μg of RBC NUDT14 (GST). MW markers (right) 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

