

Nudix (Nucleoside Diphosphate linked moiety X)-type Motif 12 (NUDT12) (His)

**CATALOG NO.:** NUC-11-342

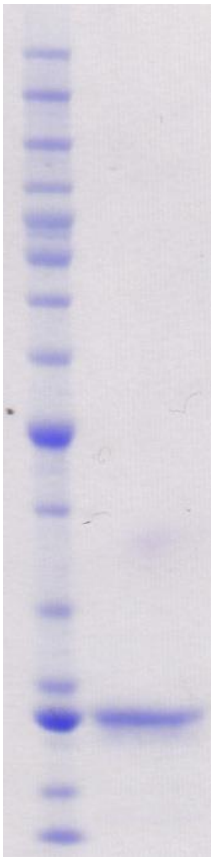
**LOT NO.:**

**DESCRIPTION:** Human recombinant NUDT12 (residues 312-462; Genbank Accession #NM\_031438.2; MW = 20.0 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, 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  $\mu\text{l}$ ) or storage of diluted enzyme is not recommended.



**Coomassie blue-stained SDS-PAGE (12% acrylamide) of 2 $\mu\text{g}$  of RBC NUDT12 (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.