

Nudix (Nucleoside Diphosphate linked moiety X)-type Motif 5 (NUDT5) (GST)

**CATALOG NO.:** NUC-11-345

**LOT NO.:**

**DESCRIPTION:** Human recombinant NUDT5 (residues 2-219; Genbank Accession #NM\_014142.2; MW = 51.1 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

**PURITY:** >85% 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 NUDT5 (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.