

Nudix (Nucleoside Diphosphate linked moiety X)-type Motif 16 (NUDT16), transcription variant 2 (GST)

**CATALOG NO.:** NUC-11-344

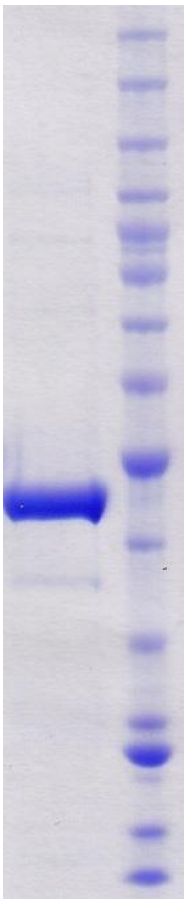
**LOT NO.:**

**DESCRIPTION:** Human recombinant NUDT16 (residues 2-195; Genbank Accession #NM\_152395.2; MW = 48.0 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  $\mu\text{l}$ ) or storage of diluted enzyme is not recommended.



**Coomassie blue-stained SDS-PAGE (12% acrylamide) of 5 $\mu\text{g}$  of RBC NUDT16 (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.