

NUDT14 (His) (UDPG pyrophosphatase)

**CATALOG NO.:** NUC-11-309

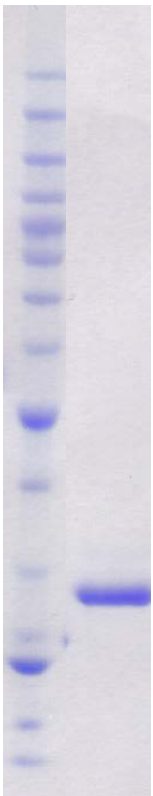
**LOT NO.:**

**DESCRIPTION:** Human recombinant NUDT14 (nudix (nucleoside diphosphate linked moiety X)-type motif 14) protein (residues 2-222; Genbank Accession #NM\_177533.4; MW = 27.0 kDa) expressed as an N-terminal His-fusion protein in *E. coli*.

**PURITY:** >90% by SDS-PAGE

**SUPPLIED AS:** µg/µL 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 4.6 µg of RBC NUDT14 (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.