

NUDT16 (His) (IDPase; U8 snoRNA-decapping enzyme)

CATALOG NO.: NUC-11-310

LOT NO.:

DESCRIPTION: Human recombinant nudix (nucleoside diphosphate linked moiety X)-type motif 16 (NUDT16) , (residues 2-195; Genbank Accession # NM_152395.2; MW = 24.2 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 4.2 µg of RBC NUDT16 (His) 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.