

OGT

O-linked N-acetylglucosamine (GlcNAc) transferase (UDP-N-acetylglucosamine:polypeptide-N-acetylglucosaminyl transferase) (OGT), transcript variant 1, (His-FLAG)

CATALOG NO.: MSC-21-176**LOT NO.:****DESCRIPTION:** Human recombinant OGT (residues 382-1046; Genbank Accession # NM_181672.2; MW = 121.2 kDa) expressed with an N-terminal His-FLAG-tag in Sf9 insect cells.**PURITY:** >95% by SDS-PAGE**SUPPLIED AS:** $_ \mu\text{g}/\mu\text{L}$ in 50 mM HEPES, pH 7.1, 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 2 μg of RBC OGT (His-FLAG). 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.

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com