

Recombinant OGT protein

Catalog No: 31524, 31924

Lot No: 19416001

Expressed In: Baculovirus

Quantity: 20 µg

Concentration: 0.55 µg/µl

Source: Mouse

Buffer Contents: Recombinant full length OGT protein is supplied at a concentration of 0.55 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

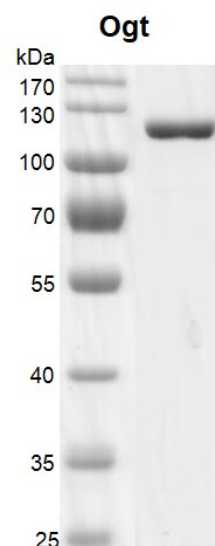
Background: OGT (O-linked N-acetylglucosamine (GlcNAc) transferase) catalyzes the transfer of a single N-acetylglucosamine from UDP-GlcNAc to serine or threonine residues in cytoplasmic and nuclear proteins resulting in their modification with a beta-linked N-acetylglucosamine (O-GlcNAc). OGT can glycosylate a large and diverse number of proteins including HCF1, AKT1, MLL5 and histone H2B. It can regulate their cellular processes via cross-talk between glycosylation and phosphorylation or by affecting proteolytic processing. Specifically OGT has been shown to interact directly with TET2 and TET3, enzymes that catalyze the oxidation of 5-methylcytosine on DNA. The TET/OGT interaction leads to GlcNAc acetylation of HCF1, a protein component of the COMPASS complex, which regulates H3K4 methylation and gene expression.

Protein Details: Recombinant mouse OGT was expressed in a baculovirus expression system as the full length protein (accession number NP_001277464.1) with an N-terminal FLAG tag. The molecular weight of the protein is 118.7 kDa. The purity of OGT protein is >90 by SDS-PAGE.

Application Notes: This protein is useful for the study of protein-protein interaction, enzyme kinetics, screening inhibitors, and selectivity profiling.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



Recombinant OGT protein gel.
OGT protein was run on an 8% SDS-PAGE gel and stained with Coomassie Blue.