

## O-GlcNAc antibody (mAb)

---

**Catalog Nos:** 61453, 61454

**RRID:** AB\_2793643

**Clone:** CTD110.6

**Isotype:** IgM

**Application(s):** IP, WB

**Reactivity:** Not Species Specific

**Quantities:** 100 µg, 10 µg

**Purification:** IgM purification

**Host:** Mouse

**Concentration:** 1 µg/µl

---

**Background:** **O-GlcNAc** ( $\beta$ -O-linked N-acetylglucosamine) is a ubiquitous and dynamic post-translational modification present on numerous cytoplasmic and nuclear proteins and is thought to play a regulatory role in numerous cellular processes including homeostatic mechanisms and gene expression. OGT (O-linked N-acetylglucosamine (GlcNAc) transferase) catalyzes the transfer of a single N-acetylglucosamine from UDP-GlcNAc to a serine or threonine residues whereas OGA / O-GlcNAcase (O-linked N-acetylglucosamine (GlcNAc)-ase) is a glycosidase that removes O-GlcNAc. O-GlcNAc modifications have been described to directly regulate the activities of a variety of transcription factors including Sp1, estrogen receptors, STAT5, NF- $\kappa$ B, p53, YY1, E1f-1, c-Myc, Rb, PDX-1, CREB, fork-head, and others. In addition, RNA polymerase II and the basal transcription factors are also extensively modified by O-GlcNAc. O-GlcNAc modification has been reported to suppress or enhance transcription, depending on the promoter and other associated proteins.

**Immunogen:** This antibody was raised against a peptide containing O-GlcNAc - serine. This antibody recognizes single O-GlcNAc residues in  $\beta$ -O-glycosidic linkage to serine and threonine.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at  $-20^{\circ}\text{C}$  for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

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