

OGA/O-GlcNAcase antibody (pAb)

Catalog Nos: 61425, 61426

RRID: AB_2793630

Isotype: IgG

Application(s): IP, WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 115 kDa

Background: **OGA / O GlcNAcase** (O-linked N-acetylglucosamine (GlcNAc) -ase) is a glycosidase that removes O-linked N-acetylglucosamine (O-GlcNAc) modifications from serine and threonine residues on nuclear and cytoplasmic proteins. This dynamic modification is catalyzed by OGT and plays a role in homeostatic mechanisms and influences gene expression.

Immunogen: This antibody was raised against a peptide within the N-terminal region of human OGA/ O-GlcNAcase.

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

Application Notes:

Applications Validated by Active Motif:

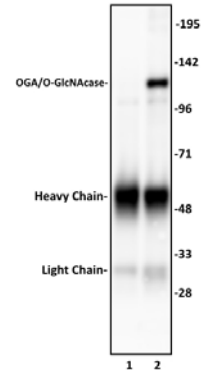
IP: 10 µl per IP

WB: 1:500 - 1:2,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

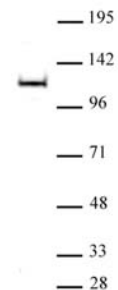
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



OGA/O-GlcNAcase antibody (pAb) tested by Immunoprecipitation.

10 µl of OGA/O-GlcNAcase antibody was used to immunoprecipitate OGA/O-GlcNAcase from 400 µg of HeLa whole cell extract (lane 2). 10 µl of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the OGA/O-GlcNAcase antibody at a dilution of 1:1,000.



OGA/O-GlcNAcase pAb tested by Western blot.

OGA/O-GlcNAcase detection by Western blot. The analysis was performed using 20 µg HeLa whole cell extract and OGA/O-GlcNAcase at a 1:1,000 dilution.