# O-GlcNAc monoclonal antibody (CTD110.6) (purified)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ENZ-ABS112A-0100

100µg

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

# Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Nameβ-O-linked N-acetylglucosamine

**Application** IF, IHC (PS)

**Application Notes**Milk proteins may cross-react with the antibody therefore

we recommend using BSA or goat serum as a blocker.

Clone CTD110.6

**Crossreactivity** Does not cross-react with peptide determinants or other

closely-related carbohydrate antigens.

**Formulation** Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

**Host** Mouse

Immunogen Synthetic peptide containing serine-O-GlcNAc.

lsotype IgM

Purity Detail Purified by affinity chromatography.

**Recommendation Dilutions/Conditions**For immunofluorescence microscopy, a concentration

range of 2.5 – 10 μg/ml is recommended. For

immunohistochemical staining on formalin-fixed paraffinembedded tissue sections, the suggested use of this reagent is  $1.0-10~\mu g/ml$ . It is recommended that the reagent be titrated for optimal performance for each

application.

Species Reactivity Species independent

### **Specificity**

Recognizes Ser-O-GlcNAc and Thr-O-GlcNAc in eukaryotes. Interaction between the antibody and the Seror Thr-O-GlcNAc epitope can be blocked by competition with free GlcNAc.

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-