

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

This antibody is covered by our [Worry-Free Guarantee](#).

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 Name β -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.

Isotype IgM

Purity Detail Purified by affinity chromatography.

Recommendation Dilutions/Conditions For immunofluorescence microscopy, a concentration range of 2.5 – 10 μ g/ml is recommended. For immunohistochemical staining on formalin-fixed paraffin-embedded tissue sections, the suggested use of this reagent is 1.0 – 10 μ 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 Ser- or 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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)