HA-tag polyclonal antibody (biotin conjugate)

Epitope tags provide a method to localize gene products in a variety of cell types, study the topology of proteins and protein complexes, identify associated proteins, and characterize newly identified, low abundance or poorly immunogenic proteins when protein specific antibodies are not available. An HA tag is frequently incorporated into recombinant proteins for a variety of purposes. An HA-tag antibody can then be used to detect the protein when doing studies with transfected cells.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS290-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application WB

Formulation Liquid. In PBS containing 1mg/ml BSA and 0.05% sodium

azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to YPYDVPDYA of the

HA tag.

Isotype IgG

Purity Detail Antigen affinity chromatography

Species Reactivity Species independent

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024

