

Anti-HA Tag

Monoclonal Antibody Against HA tag.

Catalog # ab-hatag

For research use only

Version # 05E31-SV

PRODUCT INFORMATION

Content

- 250 µl Raw Ascites Fluid containing monoclonal Anti-HA Tag

Storage

- Product is shipped at room temperature. Anti-HA Tag should be stored at -20°C. It is recommended that the total volume be divided into useable aliquots upon receipt before freezing. Avoid repeated freeze-thaw cycles.

Description

Anti-HA Tag is a monoclonal antibody that reacts specifically with the 9 amino acid HA tag (YPYDVPDYA)

Isotype: Mouse IgG1

Concentration: 5-7 mg/ml

Quality control

This lot of Anti-HA Tag has been tested by Western Blot against various HA-tagged TLR proteins.

BACKGROUND

The monoclonal antibody Anti-HA Tag was raised against the twelve amino acid peptide CYPYDVPDYASL and recognizes the influenza hemagglutinin epitope (YPYDVPDYA) which has been used extensively as a general epitope tag in expression vectors such as the pUNO/HA-TLR plasmids. The specificity of this antibody allows for unambiguous identification of the tagged protein.

APPLICATIONS

Anti-HA Tag can be used for detection of HA-tagged proteins by Western Blot or Immunoprecipitation. The utility of this antibody for other applications has not been determined.

Different assay conditions require that serial dilution of all reagents be performed to determine optimal working concentrations. Prepare working dilution immediately before use. Storage of a working concentration may result in loss performance.

Recommended working dilution for Western Blot: 1:1000

Note: When immunoblotting whole cell extracts prepared from tissue culture, Anti-HA tag may show weak background binding to serum albumin in remnant tissue culture media. Background can be eliminated without affecting HA-specific signal by rinsing tissue culture cells with PBS prior to harvest.

References

Field J. *et al.*, 1988. Purification of RAS responsive adenyl cyclase complex from *Saccharomyces cerevisiae* by use of an epitope addition method. *Mol Cell Bio*, 8:2159-2165

Kolodziej, P., Young R., 1991. Epitope tagging and protein surveillance. *Meth. Enzymol.* 194:508-519

Bennett BD, *et al.* 2000. A furin-like convertase mediates propeptide cleavage of BACE, the Alzheimer's beta-secretase. *J. Biol. Chem.* 275(48): 37712-37717

Hellwell SB, *et al.* 2001. Components of a ubiquitin ligase complex specify polyubiquitination and intracellular trafficking of the general amino acid permease. *J. Cell Biol.* 153(4): 649-662.

TECHNICAL SUPPORT

Toll free (US): 888-457-5873

Outside US: (+1) 858-457-5873

Europe: +33 562-71-69-39

E-mail: info@invivogen.com

Website: www.invivogen.com



3950 Sorrento Valley Blvd. Suite 100
San Diego, CA 92121 - USA