

**HA.11 Clone 16B12 Monoclonal Antibody, Alexa Fluor® Labeled**  
**Catalog Number: A488-101L**  
**Available Size: 0.1 mL**

**Description:** Alexa Fluor® 488 Labeled Antibody, HA.11

**Intended Use: \*\*Research Use Only (RUO)\*\***

This product is sold for laboratory research use only, not for human or in-vivo use.

**Clone:** 16B12

**Form:** Alexa Fluor® Labeled Antibody (in PBS + 5mM Azide + 0.1% BSA)

**Host:** Mouse

**IsoType:** IgG1

**[Ab]:** 1 mg/mL

**Specificity:** Monoclonal antibody HA.11 was raised against the twelve amino acid peptide, CYPYDVPDYASL. This second-generation HA antibody is an excellent substitute for the 12CA5 monoclonal antibody. HA.11 recognizes the influenza hemagglutinin epitope (YPYDVPDYA) which has been used extensively as a general epitope tag in expression vectors. The extreme specificity of the antibody allows unambiguous identification and quantitative analysis of the tagged protein. The HA.11 antibody recognizes HA epitopes located in the middle of protein sequences as well as at the N- or C-terminus. This antibody was subsequently purified via a protein-G column and conjugated to Alexa Fluor® 488. Alexa Fluor® 488 is an excellent alternative to

fluorescein (FITC) dyes.

**Uses:** This antibody is effective in immunofluorescence (IF) and FACS.

**Suggested Working Dilution:** The optimal working dilution should be determined for each specific assay condition.

- IF: 1:1,000
- FACS: 1:1,000

**Notes:** Alexa Fluor® is a trademark of Molecular Probes, Inc.

**Storage:** Store at 4°C. Freezing temperatures will denature the product.

**References:** San-Segundo PA, Roeder GS. Role for the silencing protein Dot1 in meiotic checkpoint control. *Mol Biol Cell* 11(10):3601-3615, 2000.

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

Field J, Nikawa J, Broek D, MacDonald B, Rodgers K, Wilson I, Lerner R, Wigler M. Purification of RAS responsive adenylyl cyclase complex from *Saccharomyces cerevisiae* by use of an epitope addition method. *Mol Cell Biol* 8:2159-2165, 1988.

**Warranty/Conditions:** Covance products may not be resold or modified for resale without prior written approval.

**Related Products:**

- HA.11 Monoclonal Antibody, Purified  
Catalog Number MMS-101P

- HA.11 Monoclonal Antibody  
Catalog Number MMS-101R
  
- HA.11 Polyclonal Antibody  
Catalog Number PRB-101C
  
- HA.11 Polyclonal Antibody, Purified  
Catalog Number PRB-101P

**Product Revision Date: 2/24/2009**