HA.11 monoclonal antibody (16B12)

This antibody is covered by our Worry-Free Guarantee.

Citations: 8

View Online »

Ordering Information

Order Online »

ENZ-ABS118-0500	0.5ml
ENZ-ABS118-1000	1ml
ENZ-ABS118-0200	200µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Hemagglutinin tag

Application IF, IHC, IP, WB

Clone 16B12

Formulation Liquid. Ascites.

Host Mouse

Immunogen Twelve amino acid peptide CYPYDVPDYASL.

lgG1

Specificity Recognizes the influenza hemagglutinin epitope

(YPYDVPDYA) which has been used extensively as a general epitope tag in expression vectors. Recognizes HA epitopes located in the middle of protein sequences as

well as at the N- or C-terminus.

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

.

Last modified: May 29, 2024

