## HLA-E (human) monoclonal antibody (MEM-E/02)

This antibody is covered by our Worry-Free Guarantee.

Citations: 3

View Online »

**Ordering Information** 

Order Online »

ALX-805-701-C100

100µg

Manuals, SDS & CofA

**View Online** »

## **Handling & Storage**

**Use/Stability** Stable for at least 6 months when stored at +4°C.

**Handling** Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Application IHC (PS), WB

Clone MEM-E/02

**Crossreactivity** Does not cross-react with HLA-A, -B, -C or -G.

**Formulation** Liquid. In PBS, pH 7.4, containing 15mM sodium azide.

**Host** Mouse

**Immunogen** Recombinant human HLA-E denatured heavy chain.

lgG1

Purity Detail Protein A-affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (paraffin sections, recommended

starting dilution is 5-10µg/ml; heat antigen retrieval (sodium citrate); incubate 1 hour at room temperature; detection DAB)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Species Reactivity Human

**Specificity** Recognizes denatured heavy chain of human HLA-E.

Specificity was confirmed on HLA-G/HLA-E Workshop

(Victoria 2002).

P13747

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com