

# 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-805-7010	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 6 months when stored at +4°C.
<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Application</b>	IHC (PS), WB
<b>Clone</b>	MEM-E/02
<b>Crossreactivity</b>	Does not cross-react with HLA-A, -B, -C or -G.
<b>Formulation</b>	Liquid. In PBS, pH 7.4, containing 15mM sodium azide.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant human HLA-E denatured heavy chain.
<b>Isotype</b>	IgG1
<b>Purity Detail</b>	Protein A-affinity purified.
<b>Recommendation Dilutions/Conditions</b>	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.
<b>Species Reactivity</b>	Human
<b>Specificity</b>	Recognizes denatured heavy chain of human HLA-E. Specificity was confirmed on HLA-G/HLA-E Workshop (Victoria 2002).

UniProt ID

P13747

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)