

# HLA-G (human) monoclonal antibody (MEM-G/9)

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

Citations: 6

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-805-700-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

# Product Details

Application ELISA, Flow Cytometry, ICC, IHC (FS), IP

Clone MEM-G/9

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

Host Mouse

Immunogen Recombinant human HLA-G refolded with β2-microglobulin and peptide.

Isotype IgG1

Purity Detail Purified by precipitation.

Recommendation Dilutions/Conditions ELISA (capture; in combination with Prod. No. ALX-804-482)Flow Cytometry (1-5µg/ml; JEG-3 cells could be used as positive control)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

Specificity Recognizes native HLA-G on the cell surface as well as soluble HLA-G molecule in its β2-microglobulin associated form.

UniProt ID P17693

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



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)