

# 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 $\beta$ 2-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 (1-5 $\mu$ g/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 $\beta$ 2-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)