

# HLA-G (human) monoclonal antibody (MEM-G/1)

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

Citations: 4

[View Online »](#)

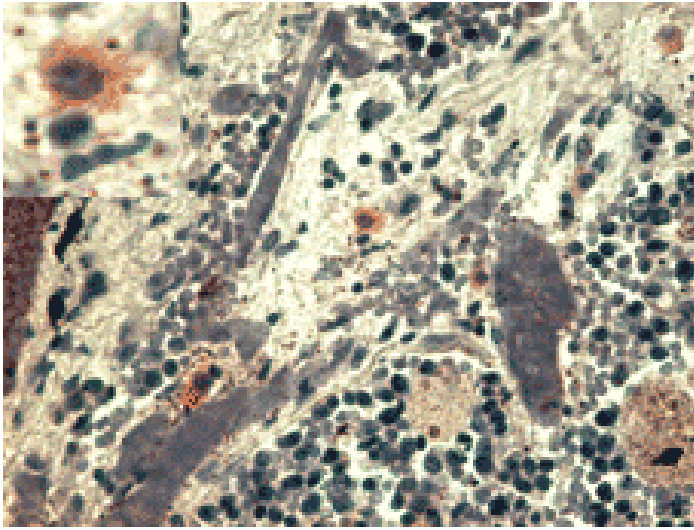
## Ordering Information

[Order Online »](#)

ALX-805-704-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Handling** Keep at +4°C. Do not freeze.

**Short Term Storage** +4°C

**Long Term Storage** +4°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** IHC (PS), WB

**Clone** MEM-G/1

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

**Host** Mouse

**Immunogen** Human recombinant HLA-G denatured heavy chain.

**Isotype** IgG1

**Purity Detail** Protein A affinity chromatography.

**Recommendation** Immunohistochemistry (paraffin sections, stain by standard ABC technique, pretraet by heat retrieval in 0.01 citrate buffer (4 x 2 minutes in microwave oven); recommended dilutions/conditions starting dilution is 1:60-1:100; incubate 1 hour at room temperature; positive tissue: Human placenta – extravillous cytotrophoblast)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 HLA-G heavy chain.

**UniProt ID** P17693



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

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

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

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

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

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

This antibody is covered by our **Worry-Free Guarantee**.