

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

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

Citations: 6

[View Online »](#)

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 β 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 μ 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 β 2-microglobulin associated form.

UniProt ID P17693

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)