

HLA-G (human) monoclonal antibody (87G)

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

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-805-719-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ELISA, Flow Cytometry, FUNC, IHC (FS)
Application Notes	Functional Application: blocks interaction of HLA-G with inhibitory receptors.
Clone	87G
Crossreactivity	Does not cross-react with mouse or rat HLA-G.
Formulation	Liquid in PBS, pH 7.4. Sterile filtered. Contains no preservatives.
Host	Mouse
Immunogen	HLA-B27 transgenic mice were immunized with H-2 identical murine cells transfected with expressing genes encoding HLA-G and human β 2-microglobulin.
Isotype	IgG2a
Purity Detail	Protein A-affinity purified.
Recommendation Dilutions/Conditions	ELISA (1 μ g/ml)Flow Cytometry (2 μ g/ml)Immunohistochemistry (frozen sections, 10 μ g/ml; incubate for 20 min. at 25°C)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human

Specificity	Recognizes membrane-bound and soluble HLA-G1 and HLA-G5.
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)