

HLA-E (human) monoclonal antibody (MEM-E/07)

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

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 6 months when stored at +4°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	Flow Cytometry, IP
Clone	MEM-E/07
Crossreactivity	Cross-reacts with classical MHC Class I molecules HLA-B7 (strongly), -B8 (moderately), -B27, -B44 (weakly).
Formulation	Liquid. In PBS, pH 7.4, containing 15mM sodium azide.
Host	Mouse
Immunogen	Recombinant human HLA-E refolded with β 2-microglobulin and peptide.
Isotype	IgG1
Purity Detail	Protein A-affinity purified.
Species Reactivity	Human
Specificity	Recognizes native surface-expressed HLA-E, but not denatured heavy chain of HLA-E.
UniProt ID	P13747
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)