

Macrophage marker monoclonal antibody (HAM56)

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|------------|----------------|
| ENZ-C34935 | ~10 µg - 0.5ml |
|------------|----------------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|--------------------|---|
| 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 | IHC (PS) |
| Clone | HAM56 |
| Crossreactivity | Cross-reacts with a subpopulation of endothelial cells including those of capillaries and smaller blood vessels. |
| Formulation | Liquid. In PBS containing 0.1% sodium azide and 1% BSA. |
| Host | Mouse |
| Immunogen | Sonicated human alveolar macrophages. |
| Isotype | IgM |
| Purity Detail | Affinity purified. |
| Source | Purified from hybridoma tissue culture supernatant. |
| Species Reactivity | Human, Monkey |
| Specificity | Recognizes fixed macrophages such as “tingible macrophages”, interdigitating macrophages of lymph nodes or tissue macrophages. |
| 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)