

# LYVE-1 (mouse), pAb

Useful to distinguish lymphatic invasion by malignant tumor cells from blood vessel invasion.

## Ordering Information

[Order Online »](#)

RLT-103-PA50AG	50µg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

Application	ELISA, Flow Cytometry, WB
Formulation	Lyophilized from PBS, pH 7.4. Contains no preservatives.
Host	Rabbit
Immunogen	Recombinant mouse soluble LYVE-1 (aa 24-228).
Isotype	IgG
Purity Detail	Antigen-affinity purified.
Reconstitution	Reconstitute with sterile water (>0.5 mg/ml).
Species Reactivity	Mouse
Technical Info / Product Notes	For the Original Manufacturer's data sheet please <a href="#">click here</a> .



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)