

LYVE-1 (mouse), pAb

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

Ordering Information

[Order Online »](#)

RLT-103-PA50	200µg
--------------	-------

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	Protein A 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 click here .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com