

# MVP/LRP (human) monoclonal antibody (LRP-56)

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

Citations: 11

[View Online »](#)

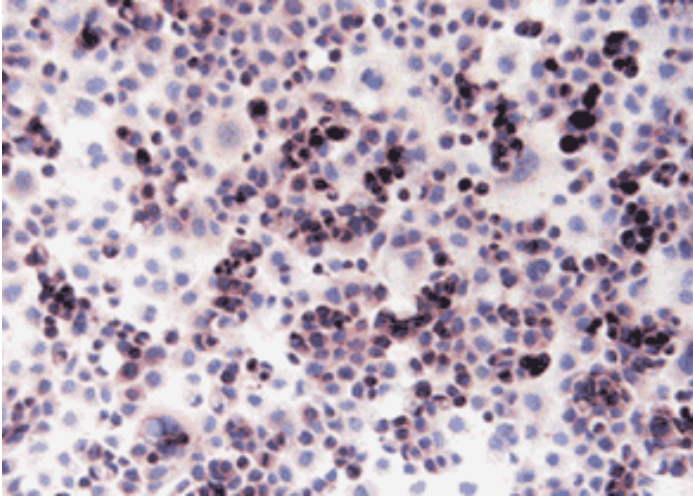
## Ordering Information

[Order Online »](#)

ALX-801-005-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



**Figure:** Liver cells stained with LRP-56 (ALX-801-005).

## Handling & Storage

**Short Term Storage** +4°C

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Major vault protein/Lung resistance protein

**Application** Flow Cytometry, ICC, IHC (FS), IHC (PS)

**Application Notes** Flow Cytometry: cell permeabilization required)  
Immunocytochemistry: cytospin preparations

**Clone** LRP-56

**Formulation** Liquid. In serum free culture supernatant containing 0.7% BSA and 0.1% sodium azide.

**Host** Mouse

**Immunogen** Human MVP/LRP (major vault protein/lung resistance protein).

**Isotype** IgG2b

**Positive Control** Cell lines: Appropriate drug-sensitive parental cell lines and their multidrug resistant derivatives prepared by the same methods as the test specimens should be used as control materials. Tissue: Human colon, kidney, bronchus or adrenal are recommended.

**Recommendation Dilutions/Conditions** Recommended starting dilution is 1:20. Suggested dilutions/conditions may not be available for all applications. Optimal should be determined individually for each application.

**Species Reactivity** Human

<b>Specificity</b>	Recognizes an internal epitope (P110) of human MVP/LRP, which is overexpressed in various human non-P-glycoprotein MDR tumor cell lines.
<b>UniProt ID</b>	Q14764
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .

**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)