

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 »](#)

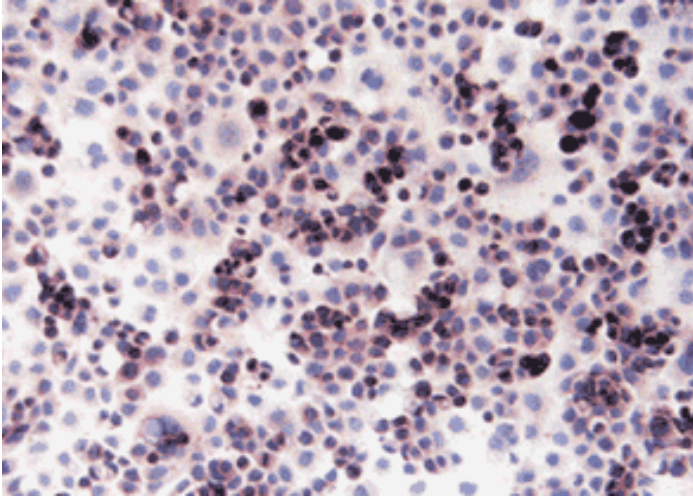


Figure: Liver cells stained with LRP-56 (ALX-801-005).

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

Specificity Recognizes an internal epitope (P110) of human MVP/LRP, which is overexpressed in various human non-P-glycoprotein MDR tumor cell lines.

UniProt ID Q14764

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)