

# Breast cancer resistance protein monoclonal antibody (BXP-53)

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

Citations: 31

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-801-036-C050	50µg
ALX-801-036-C100	100µg

Manuals, SDS & CofA

[View Online »](#)

# 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 BCRP, ABCG2, MXR, CD338

Application ICC, IHC (FS), IHC (PS), WB

Clone BXP-53

Formulation Liquid. In PBS, containing 0.1% BSA, and 0.02% sodium azide.

Host Rat

Immunogen Recombinant mouse BCRP (breast cancer resistance protein; MXR) (aa 221-394).

Isotype IgG2a

Recommendation Dilutions/Conditions Immunocytochemistry: Use 1:20-1:50 dilution on acetone fixed cytospin preparations.Immunohistochemistry (frozen sections, paraffin sections): Use 1:20 on acetone fixed frozen sections followed by incubation with biotin-labelled rabbit anti-rat IgG and HRP-labelled streptavidin., Western Blot: Use 1:20-1:50 dilution and HRP-labelled anti-rat IgG.Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse

Specificity Recognizes an internal epitope of BCRP.

UniProt ID Q7TMS5

Worry-free Guarantee

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

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)