

Breast cancer resistance protein (human) monoclonal antibody (BXP-21)

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

Ordering Information

[Order Online »](#)

ALX-801-029A-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 MXR, BCRP, ABCG2, CD338

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

Application Notes Immunocytochemistry: cytospin preparations

Clone BXP-21

Crossreactivity Does not cross-react with human P-glycoprotein, MRP1 or MRP2.

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

Host Mouse

Immunogen Recombinant human BCRP (breast cancer resistance protein; MXR) (aa 271-396).

Isotype IgG2a

Recommendation Dilutions/Conditions Immunocytochemistry: Use 1:20-1:50 dilution on acetone fixed cytospin preparations.Immunohistochemistry (frozen sections): Use 1:20 dilution on acetone fixed frozen sections followed by incubation with rabbit anti-mouse IgG and a monoclonal mouse APAAP complex.Immunohistochemistry (paraffin sections): Use on formaldehyde-fixed paraffin-embedded human tissues and tumors after pretreatment with 10mM citric acid, pH 6.0, for 20 min.Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human

Specificity Recognizes an internal epitope of human BCRP.

UniProt ID Q9UNQ0

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)