

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 .

Last modified: May 29, 2024



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