

Retinoblastoma protein monoclonal antibody (MAb245)

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Rb
Application	WB
Application Notes	Reported applications include ICC, IHC, and IP.
Clone	mAb245
Crossreactivity	Weakly cross-reacts with rat and mouse Rb.
Formulation	Liquid. In 10 mM PBS, pH 7.2, 150 mM NaCl, 0.09% sodium azide.
Host	Mouse
Isotype	IgG1
Species Reactivity	Human, Primate
Specificity	Recognizes underphosphorylated and highly phosphorylated Rb. Detects an epitope between aa 332-344 of human Rb.
UniProt ID	P06400
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