## Retinoblastoma protein monoclonal antibody (MAb245)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

BML-SA188-0100

100µg

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

lgG1

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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com