## CD200 recombinant monoclonal antibody (333)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

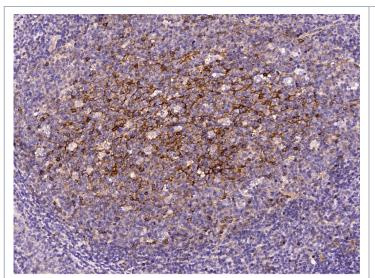
Order Online »

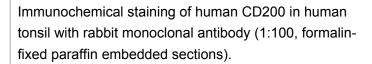
ENZ-ABS966-0100

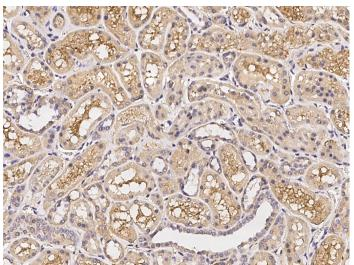
100µl

Manuals, SDS & CofA

**View Online** »







Immunochemical staining of human CD200 in human kidney with rabbit monoclonal antibody (1:100, formalinfixed paraffin embedded sections).

## **Handling & Storage**

**Handling** Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name OX-2 membrane glycoprotein

Application IHC (PS)

Clone 333

**Formulation** Liquid. 0.2 µm filtered solution in PBS. Preservative free.

**Host** Rabbit

Immunogen Recombinant human CD200 protein.

**Isotype** IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:50-

1:200)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined

individually for each application.

Species Reactivity Human

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

Last modified: September 17, 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