

CD38 recombinant monoclonal antibody (32)

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

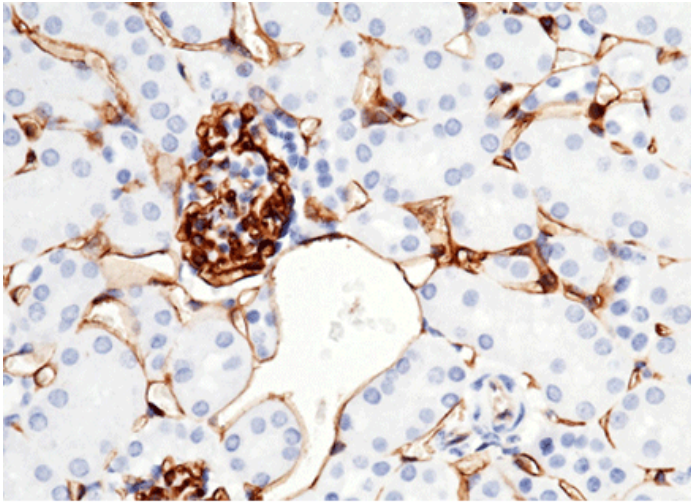
Ordering Information

[Order Online »](#)

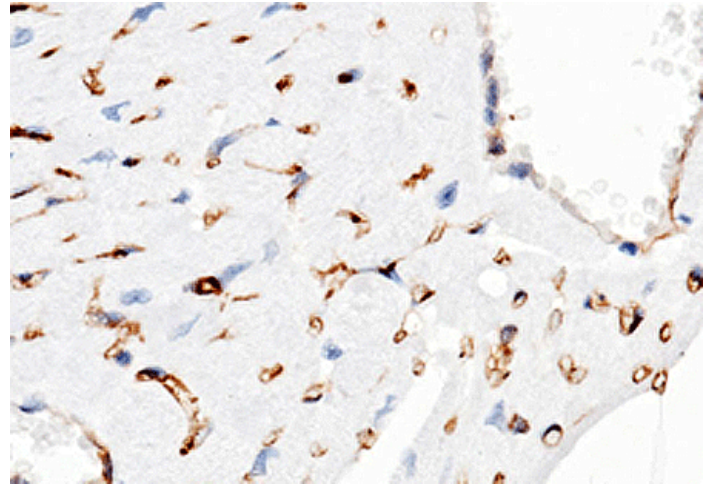
ENZ-ABS1011-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemical staining of mouse CD38 in mouse kidney with rabbit monoclonal antibody (1:500, formalin-fixed paraffin embedded sections). Positive staining was localized to endotheliocyte.



Immunohistochemical staining of mouse CD38 in mouse heart with rabbit monoclonal antibody (1:500, formalin-fixed paraffin embedded sections). Positive staining was localized to endotheliocyte.

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 ADP-ribosyl cyclase 1, Cyclic ADP-ribose hydrolase, Cyclic ADP-ribose hydrolase 1, cADPr hydrolase 1, T10, p45, NAD(+) nucleosidase

Application ELISA, IHC (PS)

Clone 32

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

Host Rabbit

Immunogen Recombinant mouse CD38 protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions ELISA (1:5'000-1:10'000)Immunohistochemistry (paraffin sections) (1:250-1:1'000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Mouse

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

.



ENZO LIFE SCIENCES,
INC.
Phone: 800 642 0430
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(EU) 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