# RAB7A polyclonal antibody

This polyclonal antibody detects RAB7A, a small GTPase involved in endosomal trafficking and autophagy.

RAB7A regulates late endosomal maturation and lysosomal biogenesis. It is essential for autophagy and degradation of cellular waste. Dysregulation of RAB7A has been linked to neurodegeneration and cancer.

This antibody is covered by our Worry-Free Guarantee.

## **Ordering Information**

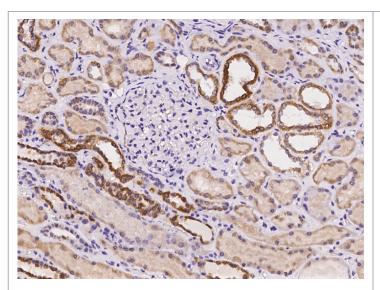
Order Online »

ENZ-ABS1071-0100

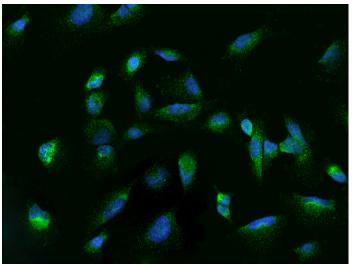
100µl

Manuals, SDS & CofA

**View Online »** 



Immunochemical staining of human RAB7A in human kidney with RAB7A polyclonal antibody at 1:100 dilution, formalin-fixed paraffin embedded sections.



Immunofluorescence staining of RAB7A in U251MG cells. Cells were fixed with 4% PFA, permeabilzed with 0.1% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit RAB7A polyclonal antibody (dilution ratio 1:100) at 4°C overnight. Then cells were stained with the Alexa Fluor®488-conjugated Goat Antirabbit IgG secondary antibody (green) and counterstained with DAPI (blue).Positive staining was localized to Cytoplasm.

### **Handling & Storage**

**Handling** Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

#### Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name Ras-related protein Rab-7a

**Application** ICC, IF, IHC (PS)

Formulation Liquid. In PBS, pH 7.0, with 0.03% Proclin300. Carrier-

free.

**Host** Rabbit

**Immunogen** E. coli-derived human RAB7A fragment.

**Isotype** IgG

Purity Detail Protein A & antigen affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (paraffin sections) (1:50-1:200)

Immunocytochemistry/Immunofluorescence (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

•



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