

# 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](#).

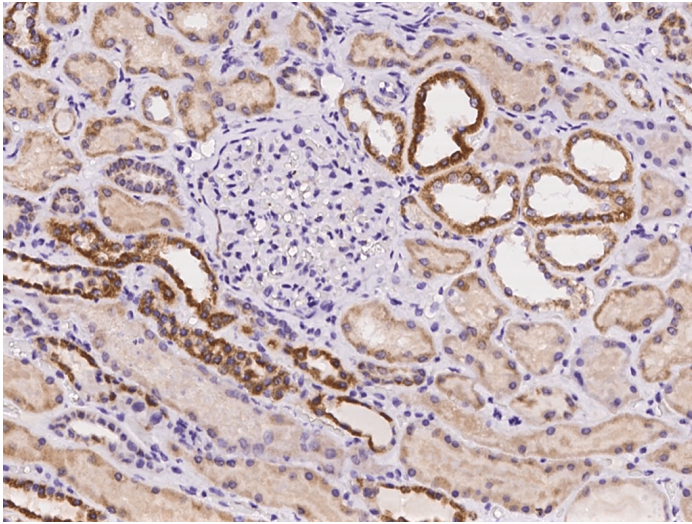
## Ordering Information

[Order Online »](#)

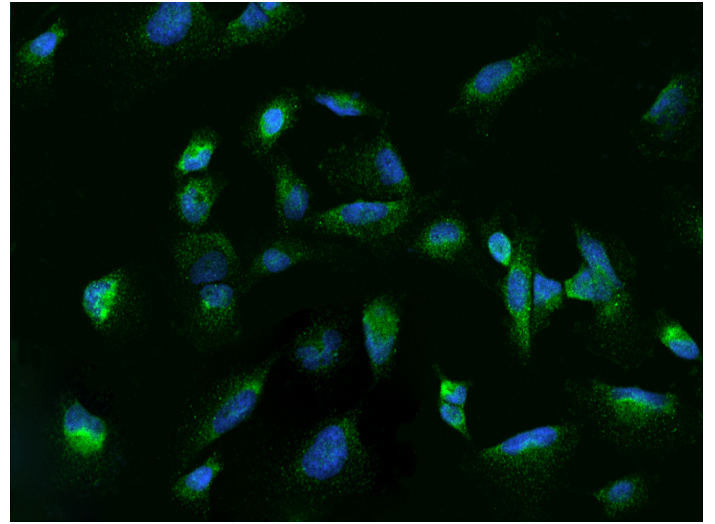
ENZ-ABS1071-0100	100µl
------------------	-------

## 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, permeabilized 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 Anti-rabbit 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	Immunohistochemistry (paraffin sections) (1:50-1:200)
Dilutions/Conditions	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 <a href="#">Worry-Free Guarantee</a> .



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)