

# Insulin polyclonal antibody

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

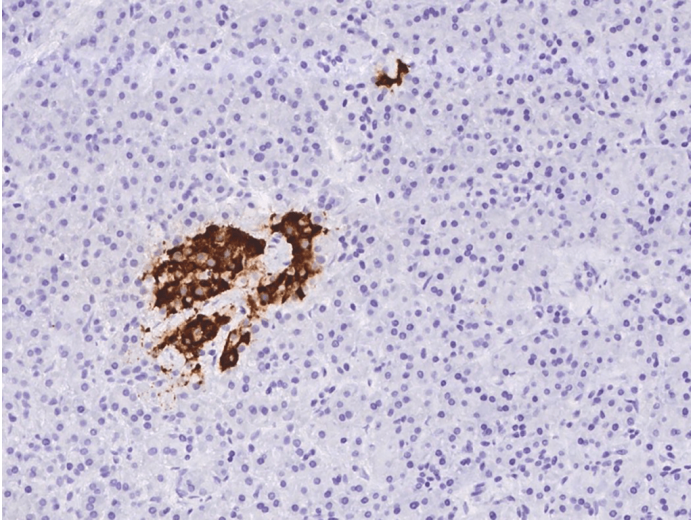
## Ordering Information

[Order Online »](#)

ENZ-ABS969-0100	100µl
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Immunochemical staining of human Insulin in human pancreas with rabbit polyclonal antibody at 1:2000 dilution, formalin-fixed 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	INS
Application	ELISA, IHC (PS)
Formulation	Liquid. 0.2 µm filtered solution in PBS. Preservative free.
Host	Rabbit
Immunogen	Recombinant human Insulin / INS protein.
Isotype	IgG
Purity Detail	Protein A and antigen affinity purified.
Recommendation Dilutions/Conditions	ELISA (1:5'000-1:10'000)Immunohistochemistry (paraffin sections) (1:500-1:5'000)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>

Last modified: September 17, 2024



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-  
fr@enzolifesciences.com](mailto:info-fr@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)