

# VISTA monoclonal antibody (027)

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

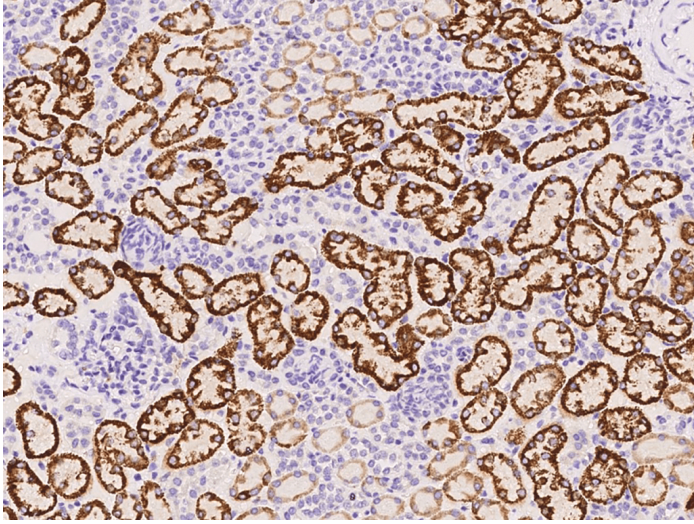
## Ordering Information

[Order Online »](#)

ENZ-ABS998-0100	100µl
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Immunochemical staining of human C10orf54 in human kidney with mouse monoclonal antibody at 1:60 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** Platelet receptor Gi24, B7-H5, V-type immunoglobulin domain-containing suppressor of T-cell activation

**Application** IHC (PS)

**Clone** 027

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

**Host** Mouse

**Immunogen** Recombinant human VISTA/GI24/B7-H5 protein.

**Isotype** IgG1

**Purity Detail** Protein A purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (paraffin sections) (1:30-1:100)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

Phone: 800.942.0430  
info-  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office

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

Belgium, The Netherlands

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

France

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

Germany

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

UK & Ireland

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