

VISTA monoclonal antibody (4C4)

VISTA/B7-H5/platelet receptor Gi24 is a single-pass type I membrane protein located at the cell surface. It is an immunoregulatory receptor which can inhibit T-cell response and may promote differentiation of embryonic stem cells, by inhibiting the BMP4 signaling pathway.

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

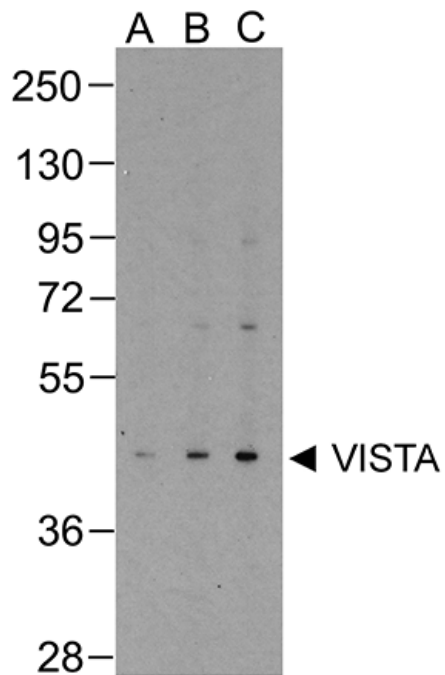
Ordering Information

[Order Online »](#)

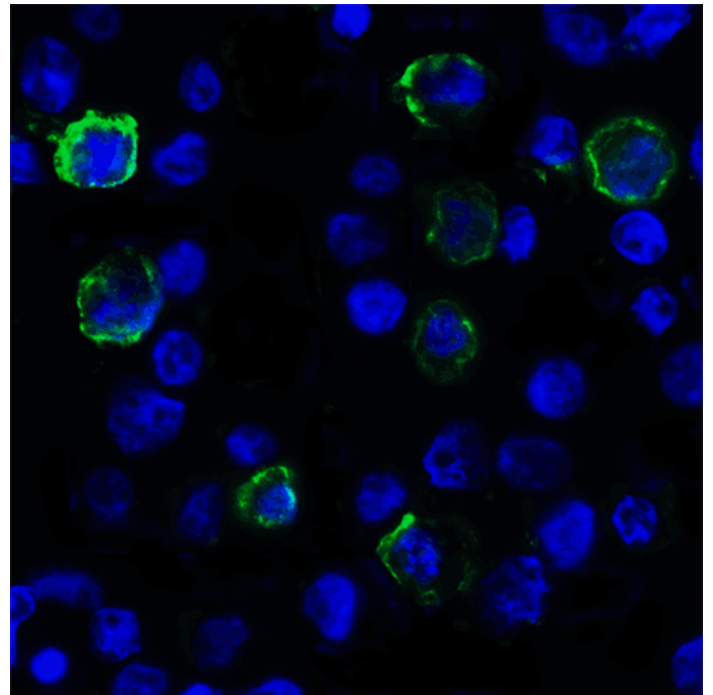
ENZ-ABS700-0100	100µg
-----------------	-------

Manuals, SDS & CofA

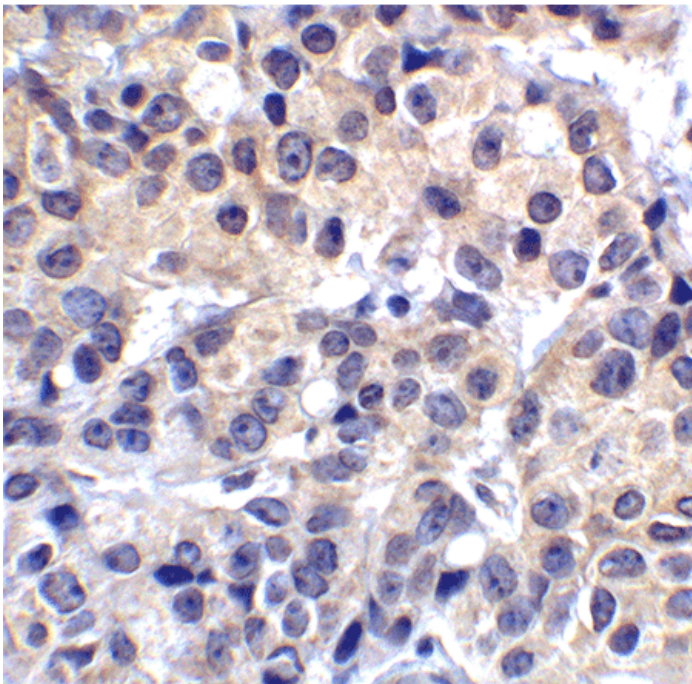
[View Online »](#)



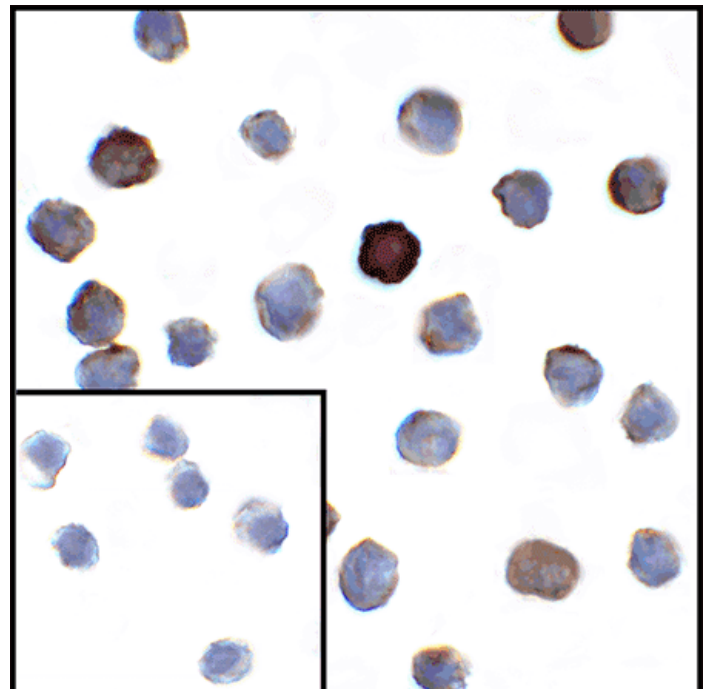
Western blot analysis of VISTA in overexpressing HEK293 cells with VISTA antibody at (A) 0.25 (B) 0.5 and (C) 1 $\mu\text{g/ml}$



Immunofluorescence of VISTA in transfected HEK293 cells with VISTA antibody at 2 $\mu\text{g/mL}$. Green: VISTA Antibody [4C4] Blue: DAPI staining



Immunohistochemistry of VISTA in human lymphoma tissue with VISTA antibody at 2 $\mu\text{g/mL}$.



Immunocytochemistry of VISTA in transfected HEK293 cells with VISTA antibody at 1 $\mu\text{g/mL}$. Lower left: Immunocytochemistry in transfected HEK293 cells with control mouse IgG antibody at 1 $\mu\text{g/mL}$.

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored as recommended.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	VSIR, V-Type immunoglobulin domain-containing suppressor of T-cell activation
Application	Flow Cytometry, ICC, IF, IHC (PS), WB
Clone	4C4
Formulation	Liquid. In PBS containing 0.02% sodium azide and 50% glycerol.
Host	Mouse
Immunogen	Extracellular domain of human VISTA.
Isotype	IgG1
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
UniProt ID	Q9H7M9
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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)