

# TIGIT monoclonal antibody (4A12)

The T cell immunoreceptor with Ig and ITIM domains (TIGIT) is a member of the PVR (poliovirus receptor) family of immunoglobulin proteins. It is expressed on several classes of T cells including follicular B helper T cells (TFH). TIGIT has been shown to bind PVR with high affinity; this binding is thought to assist interactions between TFH and dendritic cells to regulate T cell dependent B cell responses. TIGIT is upregulated on exhausted T cells in chronic viral infections and cancer. Blockade of both TIGIT and PD-1 pathways leads to tumor rejection in mice suggesting that it may be of therapeutic use against cancer.

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

Citations: 1

[View Online »](#)

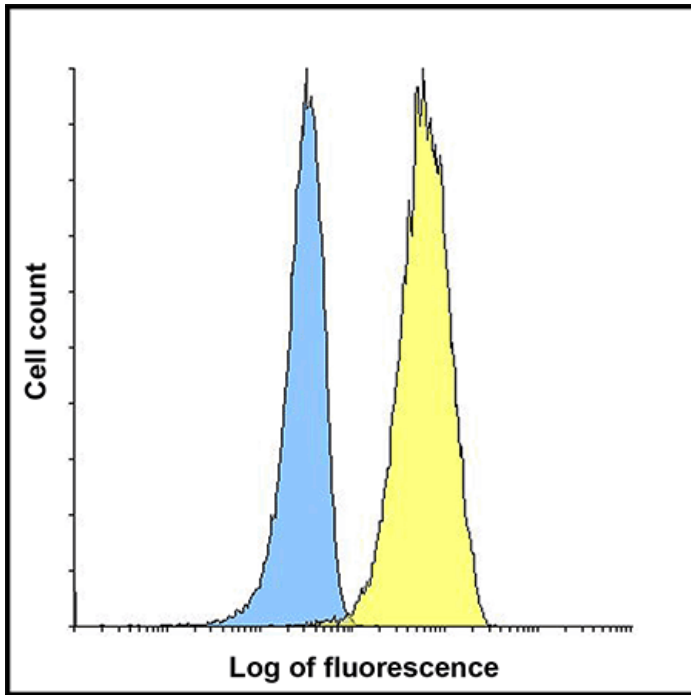
## Ordering Information

[Order Online »](#)

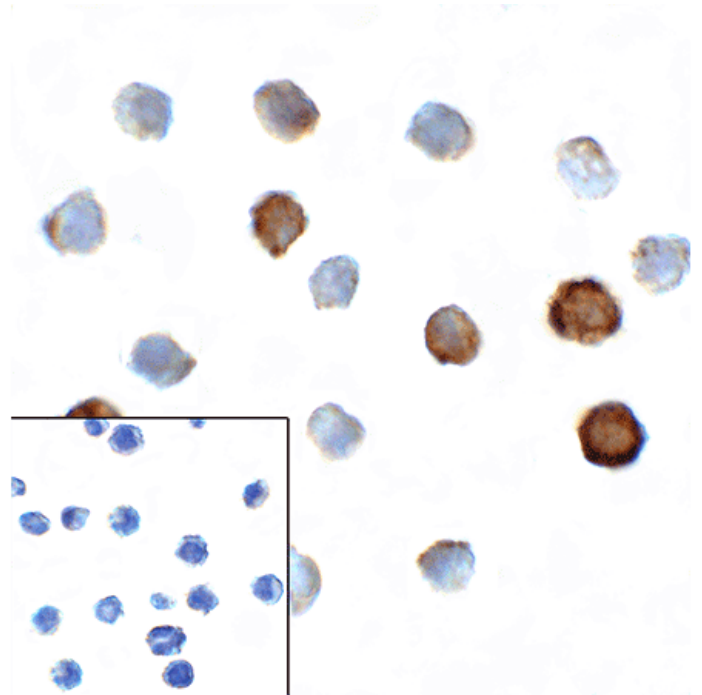
ENZ-ABS698-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

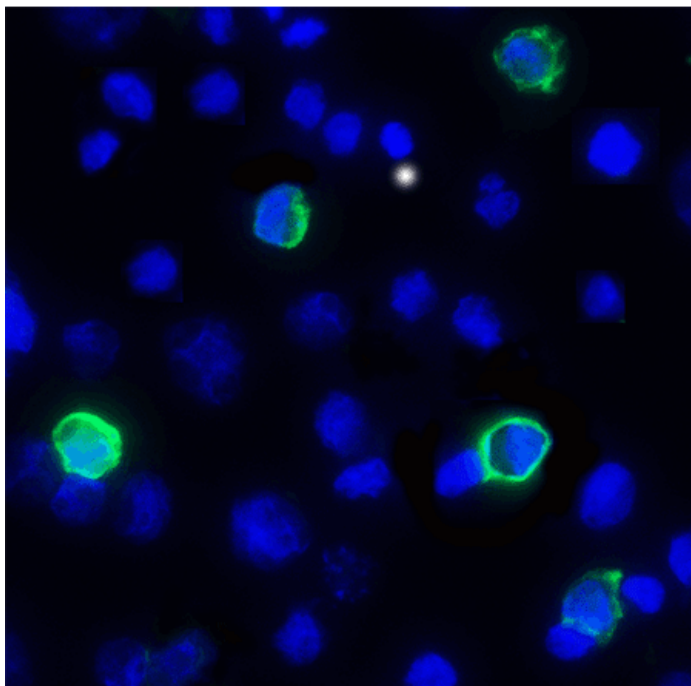
[View Online »](#)



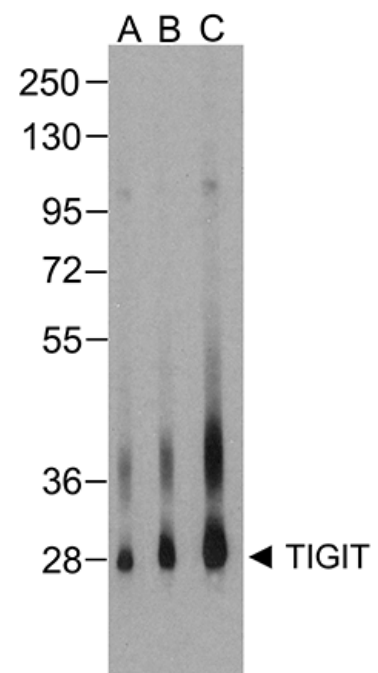
Flow cytometry analysis of TIGIT over expressing HEK293 cells using TIGIT antibody at 1  $\mu\text{g/ml}$ . Blue: untransfected HEK293 cells. Yellow: TIGIT over expressing HEK293 cells.



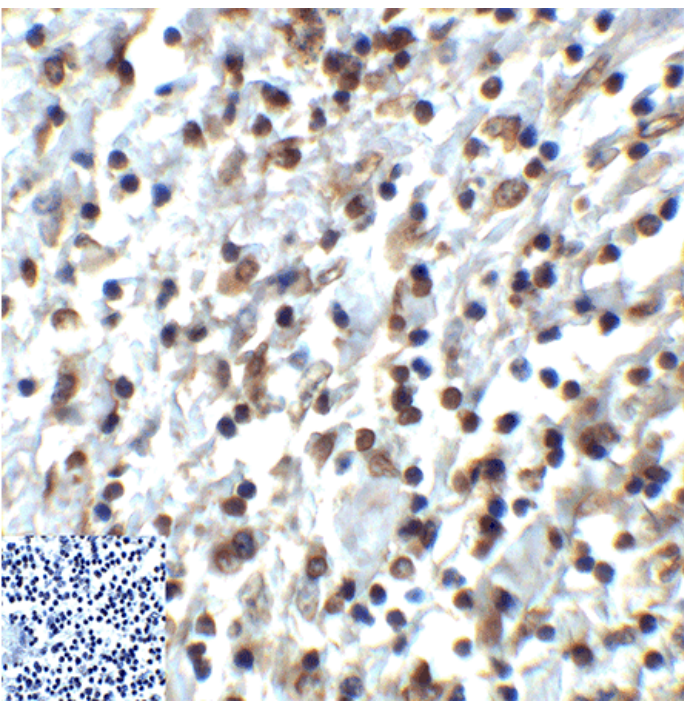
Immunocytochemistry of TIGIT in over expressing HEK293 cells using TIGIT antibody and control mouse IgG antibody (left corner box) at 1  $\mu\text{g/ml}$ .



Immunofluorescence of TIGIT in over expressing HEK293 cells using TIGIT Antibody at 1  $\mu\text{g/ml}$ . Green: TIGIT Antibody [4A12]  
Blue: DAPI staining



Western blot analysis of TIGIT in over expressing HEK293 cells using ENZ-ABS698 antibody at (A) 0.25  $\mu\text{g/ml}$ , (B) 0.5  $\mu\text{g/ml}$ , and (C) 1  $\mu\text{g/ml}$ .



Immunohistochemistry of TIGIT in human stomach carcinoma tissue using TIGIT Antibody and control mouse IgG (corner box) at 2  $\mu\text{g/ml}$ .

## 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	T-cell immunoreceptor with Ig and ITIM domains
Application	Flow Cytometry, ICC, IF, IHC (PS), WB
Clone	4A12
Formulation	Liquid. In PBS, with 0.02% sodium azide and 50% glycerol.
Host	Mouse
Immunogen	Extracellular domain of human TIGIT.
Isotype	IgG1
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
UniProt ID	Q495A1
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-  
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)