

# TMIGD2 monoclonal antibody (8A1)

TMIGD2 (transmembrane and immunoglobulin domain containing 1) is a novel adhesion molecule that is expressed in multiple tissues, primarily in cells of epithelium and endothelium origins. TMIGD2 is thought to be involved in angiogenesis and regulates cellular morphology, homophilic cell aggregation, and cell-cell interaction. TMIGD2 activity also modulates actin stress fiber formation and focal adhesion and reduces cell migration.

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

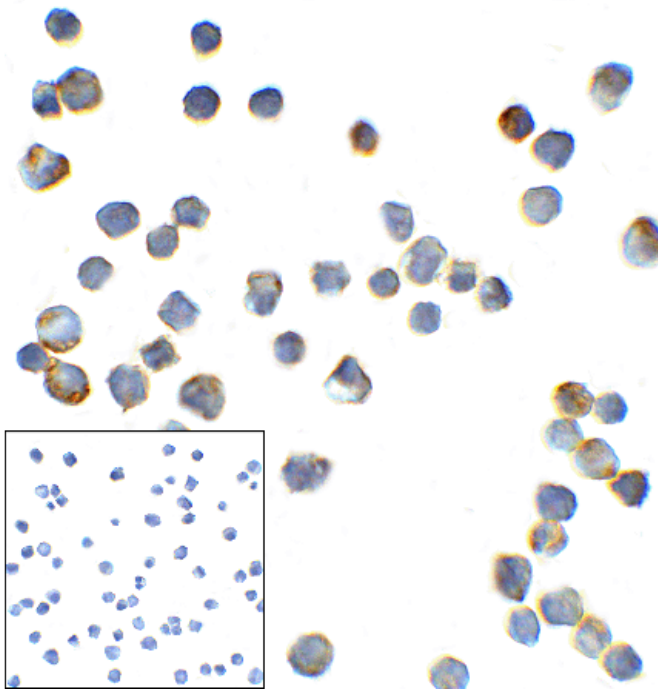
## Ordering Information

[Order Online »](#)

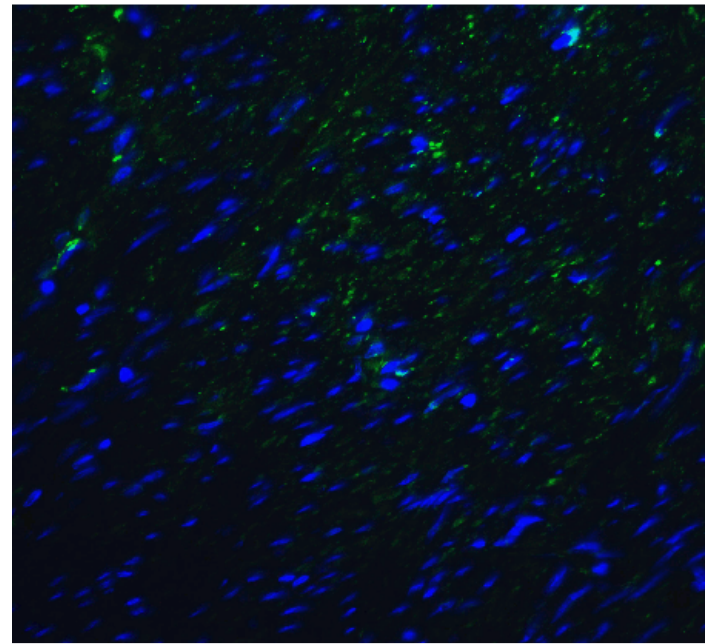
ENZ-ABS703-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

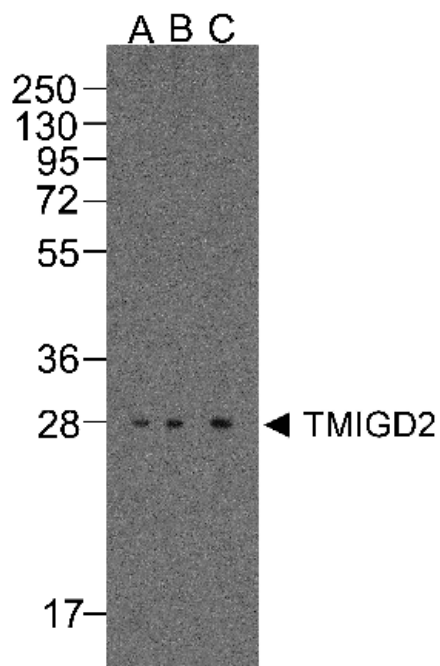
[View Online »](#)



Immunocytochemistry of TMIGD2 in transfected HEK293 cells with TMIGD2 antibody at 2  $\mu\text{g/mL}$ . Lower left: Immunocytochemistry in transfected HEK293 cells with control mouse IgG antibody at 2  $\mu\text{g/mL}$ .



Immunofluorescence of TMIGD2 in human colon carcinoma tissue with TMIGD2 antibody at 20  $\mu\text{g/mL}$ .



Western blot analysis of TMIGD2 in overexpressing HEK293 cells TMIGD2 antibody at 0.25, 0.5, and 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	IGPR-1, Transmembrane and immunoglobulin domain-containing protein 2, CD28H
Application	Flow Cytometry, ICC, IF, IHC (PS), WB
Clone	8A1
Formulation	Liquid. In PBS containing 0.02% sodium azide and 50% glycerol.
Host	Mouse
Immunogen	Extracellular domain of human TMIGD2.
Isotype	IgG1κ
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
UniProt ID	Q96BF3
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>



ENZO LIFE SCIENCES,  
Last modified **May 29, 2024**  
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-be@enzolifesciences.com](mailto:info-be@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)