

# B7-H3 monoclonal antibody (2A7)

CD276 was initially identified as a member of the B7 family of proteins. CD276 mRNA is widely expressed, but its protein expression is usually rather low. CD276 has been shown to play a role in both the costimulation as well as the coinhibition of T cell response. CD276 also plays a critical role in the control of antitumor immune responses in some cases, while in others appears to mediate antitumor immunity. It is a possible therapeutic target for at least a subset of cancers.

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

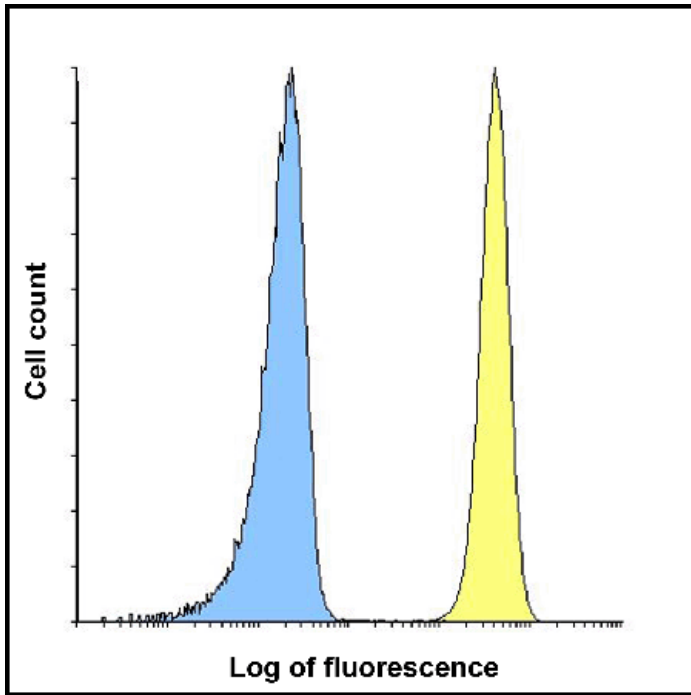
## Ordering Information

[Order Online »](#)

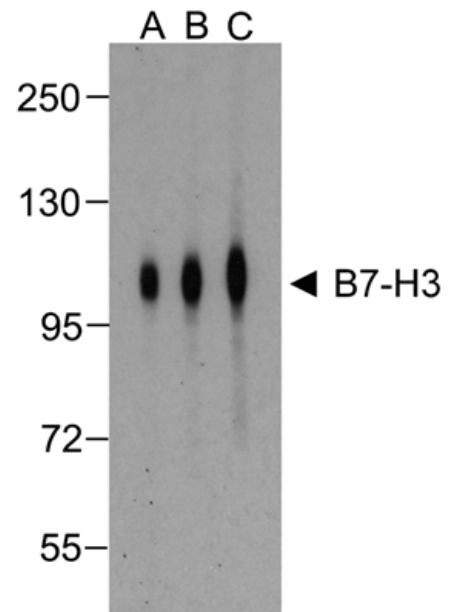
ENZ-ABS701-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

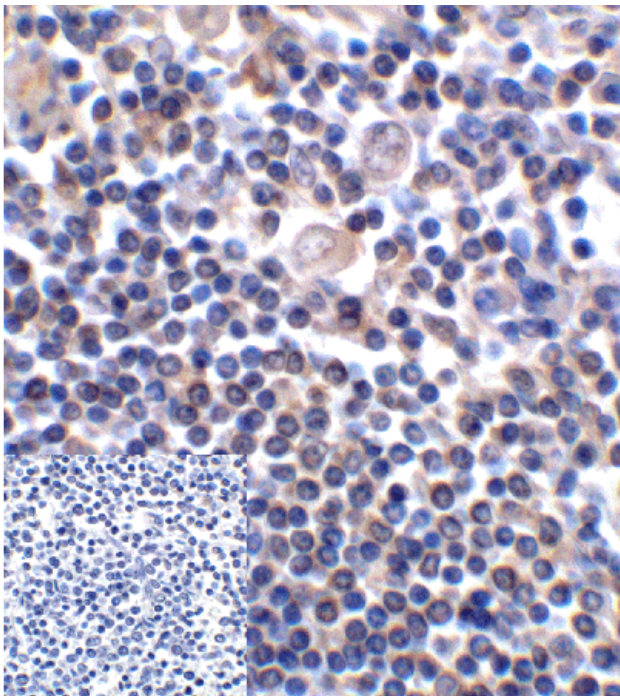
[View Online »](#)



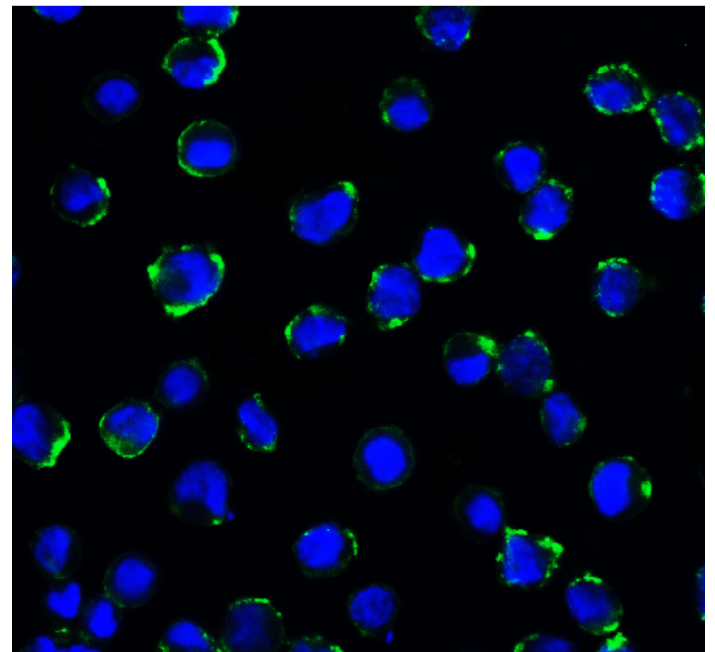
Flow cytometry analysis of B7-H3 in HEK293 cells using B7-H3 antibody at 1  $\mu\text{g/ml}$ . Blue: untransfected HEK293 cells. Yellow: B7-H3 over expressing HEK293 cells.



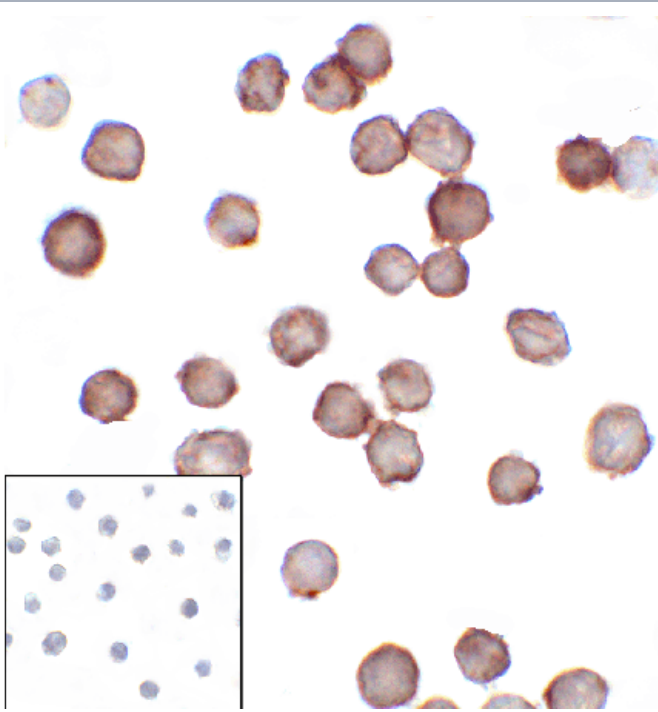
Western blot analysis of B7-H3 in HEK293 cells using B7-H3 antibody at (A) 0.25 (B) 0.5 and (C) 1  $\mu\text{g/ml}$ .



Immunohistochemistry of B7-H3 in human colon carcinoma tissue using B7-H3 Antibody and control mouse IgG (corner box) at 2  $\mu\text{g/ml}$ .



Immunofluorescence of B7-H3 in HEK293 cells using B7-H3 Antibody at 5  $\mu\text{g/ml}$ . Green: B7-H3 Antibody [2A7] Blue: DAPI staining



Immunocytochemistry of B7-H3 in HEK293 cells using B7-H3 antibody and control mouse IgG antibody (left corner box) 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	CD276
Application	Flow Cytometry, ICC, IF, IHC (PS), WB
Clone	2A7
Formulation	Liquid. In PBS containing 0.02% sodium azide and 50% glycerol.
Host	Mouse
Immunogen	Extracellular domain of human B7-H3.
Isotype	IgG1κ
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
UniProt ID	Q5ZPR3
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)