

# CD147 monoclonal antibody (BSG/7954)

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

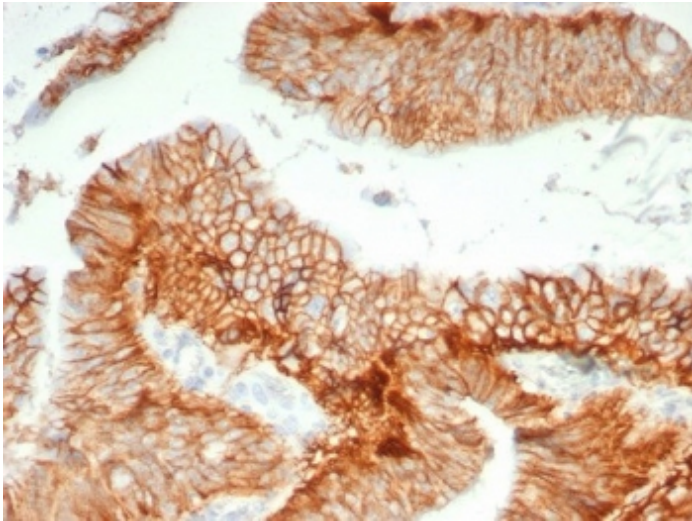
## Ordering Information

[Order Online »](#)

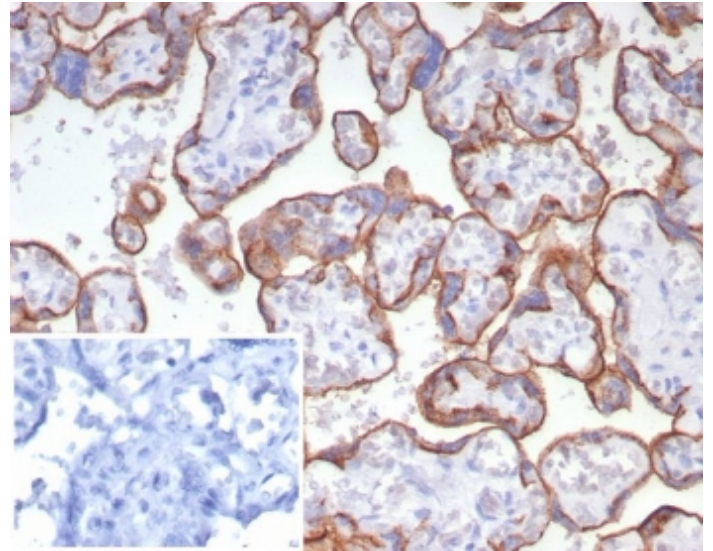
ENZ-ABS876-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

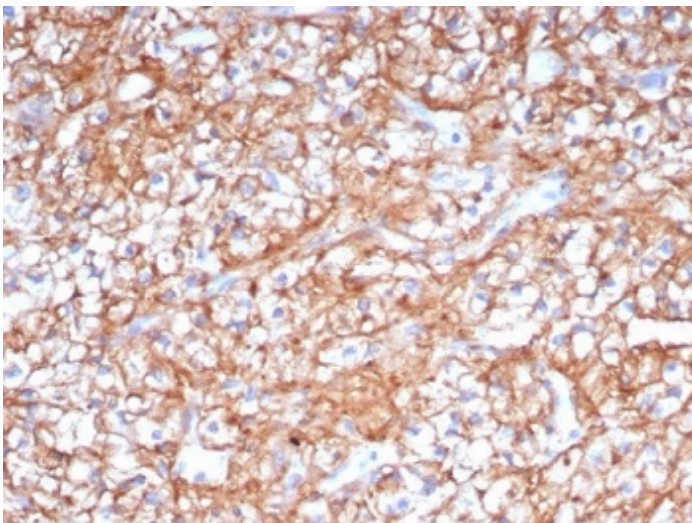
[View Online »](#)



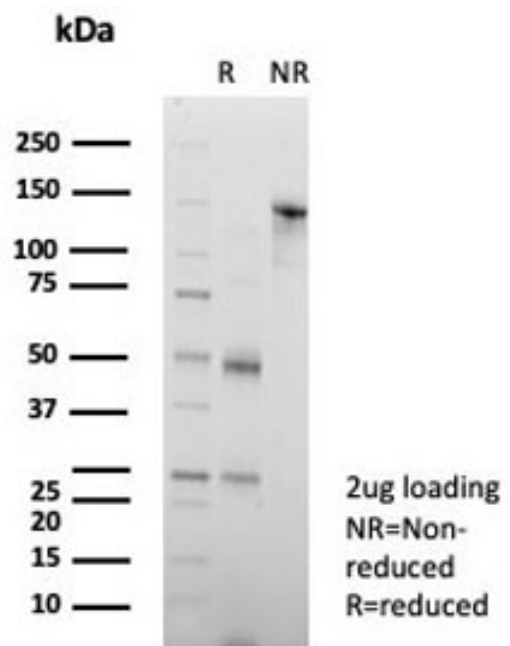
IHC staining of FFPE human colon carcinoma tissue with CD147 antibody (clone BSG/7954) at 2µg/ml. HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 20 min and allow to cool before testing.



IHC staining of FFPE human placental tissue with CD147 antibody (clone BSG/7954) at 2µg/ml. Inset: PBS used in place of primary Ab (secondary Ab negative control). HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 20 min and allow to cool before testing.



IHC staining of FFPE human renal cell carcinoma tissue with CD147 antibody (clone BSG/7954) at 2µg/ml. HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 20 min and allow to cool before testing.



SDS-PAGE analysis of purified, BSA-free CD147 antibody (clone BSG/7954) as confirmation of integrity and purity.

## Handling & Storage

Handling	Aliquot the CD147 antibody and store frozen at -20°C or colder. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	Emmprin, Basigin, Collagenase stimulatory factor
Application	ELISA, IHC (PS)
Clone	BSG/7954
Formulation	Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA (US sourced), 0.05% sodium azide.
Host	Mouse
Immunogen	Recombinant full-length human CD147.
Isotype	IgG
Purity Detail	Protein A/G affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (FFPE) (1-2µg/ml for 30 min at RT)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P35613

**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-  
be@enzolifesciences.com](mailto:info-be@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)