

# Endoglin recombinant monoclonal antibody (103)

Endoglin (CD105) is a glycoprotein expressed on the surface of endothelial cells and plays a key role in angiogenesis and cardiovascular development. It is also associated with tumor angiogenesis and is used as a marker in cancer research to study the processes of tumor growth and metastasis.

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

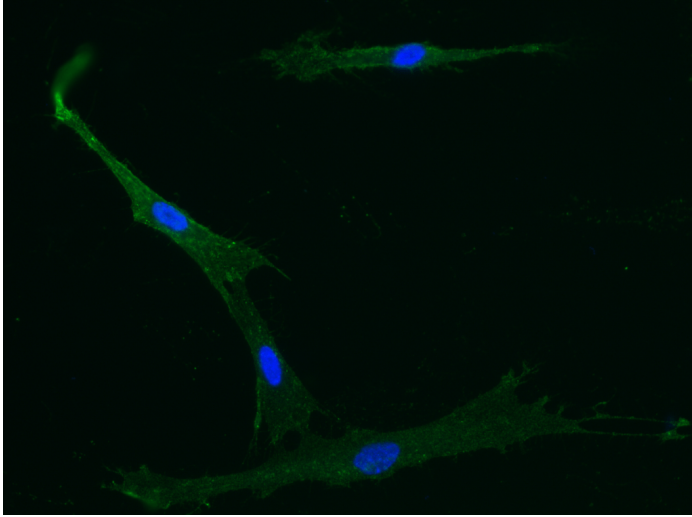
## Ordering Information

[Order Online »](#)

ENZ-ABS1030-0100	100µl
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Immunofluorescence staining of Human Endoglin/CD105 in MSC cells. Cells were fixed with 4% PFA, blocked with 10% serum, and incubated with rabbit anti-Human Endoglin monoclonal antibody (1:60) at 37°C 1 hour. Then cells were stained with the Alexa Fluor® 488-conjugated goat Anti-rabbit IgG secondary antibody (green) and counterstained with DAPI (blue). Positive staining was localized to cell membrane.

## Handling & Storage

**Handling** Avoid freeze-thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** CD105

**Application** Flow Cytometry, ICC, IF

**Clone** 103

**Formulation** Liquid. 0.2 µm filtered solution in PBS. Preservative free. Carrier-free.

**Host** Rabbit

**Immunogen** Recombinant human Endoglin / CD105.

**Isotype** IgG

**Purity Detail** Protein A purified.

**Recommendation** Immunocytochemistry / Immunofluorescence (1:20-1:100) Flow Cytometry (1:25-1:100) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

**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)