

# Thy-1 (mouse):Fc (human), (recombinant)

Thy-1 is a glycosyl phosphatidylinositol (GPI)-anchored glycoprotein of the immunoglobulin superfamily expressed in various cell types, particularly in cells of the neuronal system, activated microvascular endothelial cells, fibroblasts and blood stem cells. Thy-1 is involved in cell adhesion. Thy-1 interacts with the integrin  $\alpha$ v $\beta$ 3 (CD51/CD61) and Mac-1 (CD11b/CD18). In astrocytes, binding of Thy-1 to  $\beta$ 3-integrin promotes focal adhesion formation.

## Ordering Information

[Order Online »](#)

ALX-522-096-C050	50 $\mu$ g
------------------	------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 6 months after receipt when stored at -20°C.
<b>Handling</b>	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	CD90
<b>Concentration</b>	1mg/ml after reconstitution.
<b>Formulation</b>	Lyophilized. Contains PBS.
<b>MW</b>	~48kDa (SDS-PAGE)
<b>Purity</b>	≥90% (SDS-PAGE)
<b>Reconstitution</b>	Reconstitute with 50µl sterile distilled water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
<b>Source</b>	Produced in HEK 293 cells. The extracellular domain of mouse Thy-1 (aa 20-131) is fused to the Fc portion of human IgG1.
<b>UniProt ID</b>	P01831

Last modified: May 29, 2024



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)