

Control:Fc fusion protein (human), (recombinant)

Negative control for Fc chimeric proteins

Negative control for Fc chimeric proteins. Does not bind to a series of ligands in an ELISA assay. Does not inhibit binding between the thymoma cell line EL-4 expressing $\beta 3$ -integrin and astrocytes expressing high level of Thy-1 in mice.

Thy-1 is a glycosyl phosphatidylinositol (GPI)- anchored glycoprotein of the immunoglobulin superfamily expressed in various cell types, particularly in T-lymphocytes and cells of the neuronal system. Thy-1 is involved in cell adhesion, and the presence of a RLD motifs, which is recognized by integrins, strengthen the hypothesis that the Thy-1:integrin interaction is responsible for cell attachment. In astrocytes, binding of Thy-1 to $\beta 3$ -integrin triggers tyrosine phosphorylation of focal adhesion proteins.

Citations: 11

[View Online »](#)

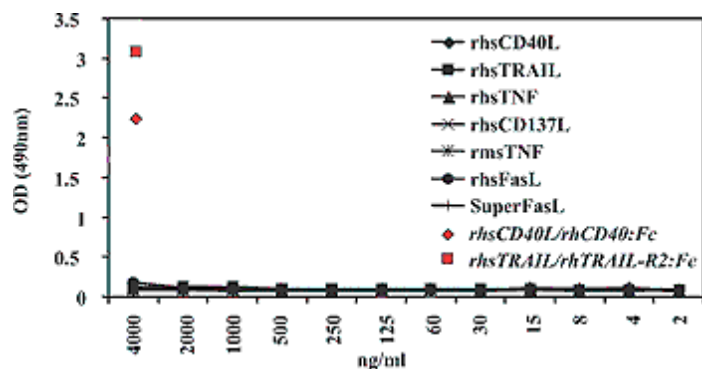
Ordering Information

[Order Online »](#)

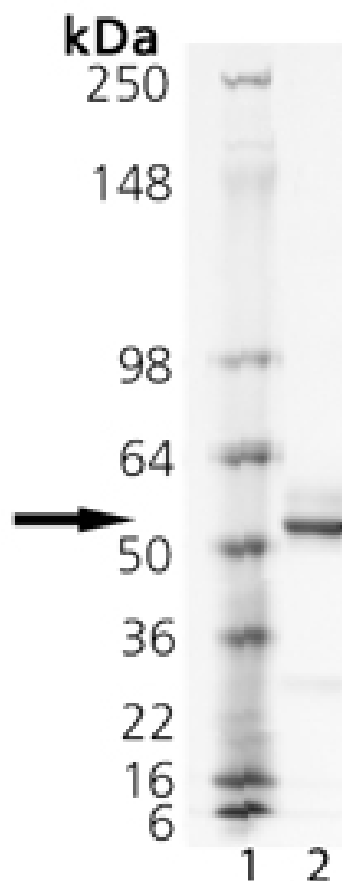
ALX-203-004-C050	50 μ g
------------------	------------

Manuals, SDS & CofA

[View Online »](#)



ELISA analysis of Control:Fc fusion protein (human), (recombinant), (Prod. No. ALX-203-004): Negative binding to members in TNF-superfamily, including, but not limited to, CD40L, TRAIL, TNF, CD137L, FasL at various concentrations as indicated. (The CD40L and TRAIL used in the assay are fully capable to bind to their native biological receptors , i.e. CD40 and TRAIL-R2, respectively, as shown by the red square and diamond.)



SDS-PAGE Analysis. Lane 1: MW Marker, Lane 2: 1 µg Fc:control protein

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored as recommended.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Thy-1 (RLE mutant)
Application Notes	Control compound for Fc Tagged Proteins, ELISA, Flow Cytometry, Functional Application (serves as a negative control in functional assays of TNF-superfamily members).
Formulation	Lyophilized. Contains PBS.
MW	~ 53 kDa
Purity	≥90% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Reconstitution	Reconstitute with 50µl sterile distilled water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
Source	Produced in HEK 293 cells. The extracellular domain of mutated hThy-1, hThy-1 (RLE) (aa 20-130), is fused to the Fc portion of human IgG1. The integrin binding sequence RLD of hThy-1 has been changed to RLE abolishing the binding of integrins.
Specificity	Does not bind to β3-integrin nor to any of a large number of tested TNF-superfamily members including rhsTRAIL, rhsCD40L, rhsAPRIL, etc.



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-
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)

Last modified: May 29, 2024

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.