

CD56 (human), (recombinant) (Fc-tag)

CD56, also known as NCAM-1, is a Protein Coding gene. CD56 is a neural adhesion protein (NCAM) that belongs to the immunoglobulin superfamily. The encoded protein is involved in cell-to-cell interactions as well as cell-matrix interactions during development and differentiation. It is involved in neuron-neuron adhesion, neurite fasciculation and the outgrowth of neurites. It has also been shown to be involved in the expansion of T cells and dendritic cells which play an important role in immune surveillance. Diseases associated with CD56 include Rabies and Bile Duct Cancer.

Ordering Information

[Order Online »](#)

ENZ-PRT325-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

kDa

M

116

66.2

45.0

35.0

25.0

18.4

14.4



As a result of glycosylation, the recombinant protein migrates as an approximately 63.8 kDa protein in SDS-PAGE under reducing conditions.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NCAM-1, Neural cell adhesion molecule 1

Endotoxin Content <1 EU/μg (LAL method)

Formulation Lyophilized from sterile 20 mM MES, 150 mM NaCl, 8% Glycerol, 10% Trehalose, pH 6.0.

MW ~83.4 kDa (SDS-PAGE)

Purity ≥95% (SDS-PAGE)

Reconstitution Reconstitute with sterile deionized water. Reconstitution instructions are lot specific.

Source Produced in HEK293 cells. A DNA sequence encoding the human CD56 (Met1-His340) was expressed with the Fc region of human IgG1 at the C-terminus.

UniProt ID P43629



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com