

CD27 (human):Fc (human), (recombinant)

The CD27 receptor (CD27, Tp55) is a member of the tumor necrosis factor receptor (TNF-R) family. CD27 is a type I transmembrane receptor, it contains 3 extracellular cysteine-rich motifs. CD27 is mainly expressed on T and B lymphocytes, and its expression is increased on T cells by the engagement of the T cell receptor complex. Engagement of CD27 induces recruitment of TNF receptor-associated factor 2 (TRAF2) and TRAF3. CD27-CD70 have been reported to be involved in several different immune mechanisms such as T, NK and B lymphocytes interactions and costimulation, early thymic differentiation, autoimmune and viral diseases and apoptosis.

Ordering Information

[Order Online »](#)

ALX-522-029-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	TNFRSF 7, CD27L receptor, T-cell activation antigen CD27
Application Notes	ELISA: binds to CD27L at 1-10 ng/ml. Can be used together with anti-human IgG1 PAb to detect membrane-bound human and mouse CD70.
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/µg purified protein (LAL test; Associates of Cape Cod).
Formulation	Lyophilized. Contains PBS.
MW	~55kDa (SDS-PAGE).
Purity Detail	Purified by single-step affinity chromatography.
Reconstitution	Reconstitute with 50µl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
Source	Produced in HEK 293 cells. The cysteine-rich region of human CD27 (aa 1-191) is fused to the Fc portion of human IgG1.
Specificity	Binds human and mouse CD70 (CD27L).
UniProt ID	P26842



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)