

CD30 (human):Fc (human), (recombinant)

The CD30 receptor (CD30, Ki-1) is a member of the tumor necrosis factor receptor (TNF-R) family. CD30 is a type I transmembrane receptor, it contains 6 extracellular cysteine-rich motifs, is shed by the TNF α -converting enzyme and is associated with the Reed-Sternberg cells of Hodgkins disease as well as with T cells and some lymphomas. Activated CD30 recruits TNF receptor-associated factor 1, 2, 3 and 5. CD30-CD30L have been reported to be involved in costimulation of T cell proliferation, induction of HIV expression, negative selection and apoptosis as well as in diabetes and several allergic, viral and autoimmune diseases.

Ordering Information

[Order Online »](#)

ALX-522-027-C050	50 μ g
------------------	------------

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	Ki-1, TNFRSF 8
Application Notes	ELISA: binds to CD30L. Can be used together with anti-human IgG1 PAb to detect membrane-bound human CD30L.
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/µg purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS.
MW	~110 kDa observed (SDS-PAGE)
Purity	≥95% (SDS-PAGE)
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 CD30 (Ki-1) (aa 1-380) is fused to the Fc portion of human IgG1 .
Specificity	Binds human CD30L (CD153).
UniProt ID	P28908



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)