

DR3 (human):Fc (human), (recombinant)

The DR3 receptor (WSL-1, Apo-3, LARD, TRAMP, TR3) is a member of the tumor necrosis factor receptor (TNF-R) family. DR3 is a type I transmembrane receptor. It is 339 aa long and contains 4 extracellular cysteine-rich motifs in its extracellular domain, while its cytoplasmic region contains a death domain. DR3 is abundantly expressed on thymocytes and T and B lymphocytes. Overexpression of DR3 mediates activation of nuclear factor NF-κB and induces apoptosis. Besides the induction of apoptosis, DR3 has been reported to regulate negative selection during thymocyte development.

Ordering Information

[Order Online »](#)

ALX-522-040-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	WSL-1, Apo-3, LARD, TRAMP, TR3, TNFRSF 12, TNFRSF 25, Death receptor 3
Application Notes	Can be used together with anti-human IgG1 PAb to detect TL1A.
Concentration	1mg/ml after reconstitution.
Formulation	Lyophilized. Contains PBS.
MW	~55kDa (under reducing conditions).
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 DR3 (aa 25-199) is fused to the Fc portion of human IgG1.
Specificity	Binds human TL1A.
UniProt ID	Q93038



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

Last modified: May 29, 2024