

# CD137 (human):Fc (human), (recombinant)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-031-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	4-1BB, TNFRSF 9
Application Notes	<p>ELISA: binds to CD137L at 10-100 ng/ml.</p> <p>Can be used together with anti-human IgG1 PAb to detect membrane-bound human CD137.</p> <p>Functional Application: Induces proliferation of monocytes.</p> <p><b>Note:</b> Induction of monocyte proliferation by CD137 required immobilization of the CD137 protein, e.g. by coating of the protein to the surface of the tissue culture dishes.</p>
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/µg purified protein (LAL test; Associates of Cape Cod).
Formulation	Lyophilized. Contains PBS.
MW	~50kDa (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 CD137 (4-1BB) (aa 1-186) is fused to the Fc portion of human IgG1.

Specificity	Binds human CD137L (4-1BBL).
UniProt ID	Q07011

Last modified: May 29, 2024



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)