

CD134L (soluble) (human), (recombinant)

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-023-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -80°C.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	OX40L, TNFSF 4, CD252
Application Notes	ELISA: binds to CD134 receptor). Can be used together with anti-FLAG [®] MAb to detect membrane-bound human CD134 (OX40).
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS.
MW	~30kDa (SDS-PAGE).
Purity Detail	Purified by single-step affinity chromatography.
Reconstitution	Reconstitute with 100μl 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 extracellular domain of human CD134L (OX40L) (aa 52-183) is fused at the N-terminus to a linker peptide (8 aa) and a FLAG [®] -tag.
Specificity	Binds to human CD134 (OX40).
Technical Info / Product Notes	Historical data indicates that CD134L activates T cell proliferation and delays neutrophil apoptosis. FLAG is a registered trademark of Sigma-Aldrich Co.

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)