

MSD[®] Human FasL Kit

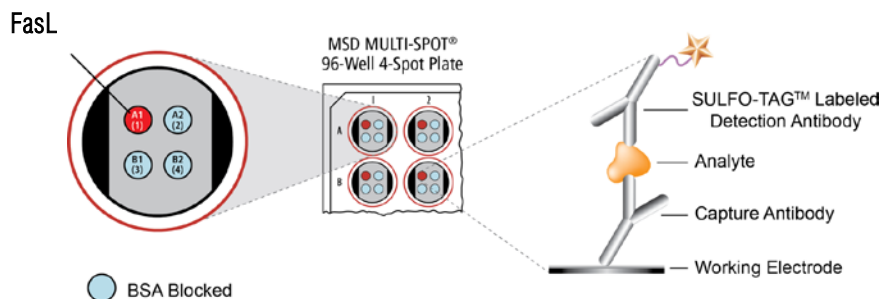
For quantitative determination in human serum, plasma, and tissue culture supernatants



Alzheimer's Disease
BioProcess
Cardiac
Cell Signaling
Clinical Immunology
Cytokines
Growth Factors
Hypoxia
Immunogenicity
Inflammation
Metabolic
Oncology
Toxicology
Vascular

Catalog Numbers

Human FasL Kit	
Kit Size	
1 plate	K151KDD-1
5 plates	K151KDD-2
25 plates	K151KDD-4



Fas Ligand (FasL or CD95L) is a 40 kDa, homotrimeric, type II membrane protein that belongs to the tumor necrosis factor family (TNF).¹ FasL is an essential effector in immune system functions; the protein induces apoptosis in Fas(CD95/APO-1)-expressing cells. In hematopoietic cells, intracellular FasL is stored in secretory lysosomes and is mobilized to the immunological synapses upon activation, a process mediated by trafficking and cytosolic reorganization adaptor proteins. FasL surface expression is regulated by post-translational ectodomain shedding by disintegrin and subsequent regulated intramembrane proteolysis, resulting in the release of soluble FasL into serum.^{2,3} FasL binding to the FasL receptor results in receptor trimerization, internalization of the Fas complex, initiation of the apoptotic signaling cascade, and cell death.⁴ FasL deficiencies have been implicated in tumor drug resistance, oncogenesis, and autoimmune disorders.⁵

The MSD Human FasL assay is available on 96-well, 4-spot plates. This datasheet outlines the performance of the assay.

Assay Sensitivity

	FasL
LLOD (pg/mL)	0.28

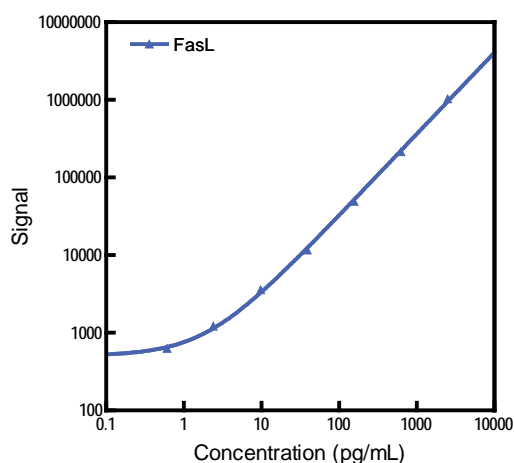
The lower limit of detection (LLOD) is a calculated concentration based on a signal 2.5 standard deviations above the background (zero calibrator blank).

Ordering information

MSD Customer Service
Phone: 1-301-947-2085
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com

Typical Standard Curve

The following standard curve is an example of the wide dynamic range of the Human FasL assay.



Conc. (pg/mL)	FasL	
	Average Signal	%CV
0	430	6.6
0.61	628	7.0
2.4	1219	3.8
9.8	3578	1.3
39	11 588	2.6
156	49 385	0.8
625	214 525	1.4
2500	1 030 610	4.2

Company Address

MESO SCALE DISCOVERY[®]
division of
Meso Scale Diagnostics, LLC.
1601 Research Boulevard
Rockville, MD 20850-3173 USA

www.mesoscale.com[®]

For Research Use Only.
Not for use in diagnostic
procedures.

MSD Biomarker Assays

Tested Samples

Normal human serum samples were diluted 2-fold and tested with the Human FasL Kit. Median and range of concentrations are displayed below. Concentrations are corrected for sample dilution.

Sample Type	Statistic	FasL
Serum	Median (pg/mL)	15.0
	Range (pg/mL)	6.4–46
	Number of Samples	20
	Samples in Quantitative Range	20

References

1. Ashkenazi A. Targeting death and decoy receptors of the tumor-necrosis factor super family. *Nat Rev Cancer*. 2002 Jun;2(6):420-430.
2. Jodo S, et al. Bioactivities of Fas Ligand-Expressing Retroviral Particles. *J Immunol* 2000 May 15;164(10):5062-5069.
3. Voss M, et al. Posttranslational regulation of Fas ligand function. *Cell Commun Signal* 2008 Dec 29;6:11.
4. Voss M, et al. Identification of SH3 domain interaction partners of human FasL (CD178) by phage display screening. *BMC Immunol* 2009 Oct 6;10:53.
5. Pitti, RM, et al. Genomic amplification of a decoy receptor for Fas ligand in lung and colon cancer. *Nature*. 1998 Dec 17;396(6712):699-703.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, WWW.MESOSCALE.COM, MSD, MSD (DESIGN), DISCOVERY WORKBENCH, QUICKPLEX, MULTI-ARRAY, MULTI-SPOT, SULFO-TAG, STREPTAVIDIN GOLD, SECTOR, SECTOR HTS, SECTOR PR, 4-SPOT (DESIGN) and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

© 2012 Meso Scale Diagnostics, LLC. All rights reserved.

For Research Use Only. Not for use in diagnostic procedures.

