

MSD[®] Human sFas Kit

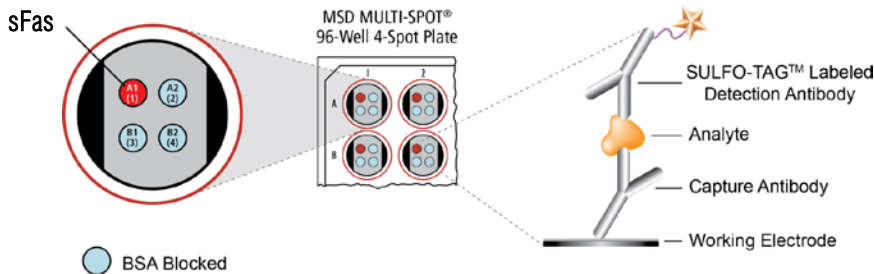
For quantitative determination in human serum, plasma, and tissue culture supernatants



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

Catalog Numbers

Human sFas Kit	
Kit Size	
1 plate	K151KBD-1
5 plates	K151KBD-2
25 plates	K151KBD-4



Soluble Fas (sFas or sCD95) is a 48 kDa, type I membrane protein that plays a critical role in apoptosis.¹ Several sFas isoforms are generated by alternative mRNA splicing; the predominant form lacks the transmembrane domain, a result of exon 6 deletion, while less active isoforms lack combinations of exons 3, 4, 6, and 7. Functionally, sFas inhibits extracellular Fas/FasL binding, impairing the homeostatic regulation of immune responses such as Fas-induced apoptosis.²

sFas is present in activated human lymphocytes and tumor cell lines supernatants. Elevated levels have been observed in patients suffering from systemic lupus erythematosus and hematopoietic malignancies, suggesting that sFas modulates autoimmune diseases and oncogenesis.³⁻⁴

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

Assay Sensitivity

	sFas
LLOD (pg/mL)	0.44

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

Company Address

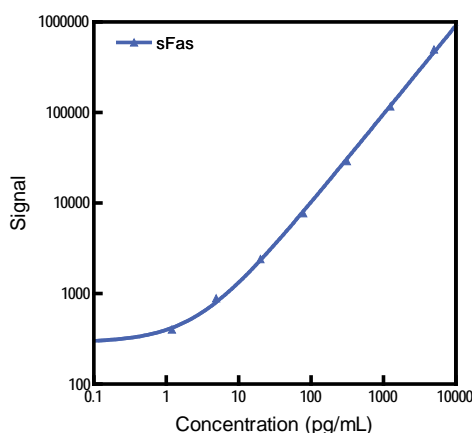
MESO SCALE DISCOVERY[®]
division of
Meso Scale Diagnostics, LLC.
1601 Research Boulevard
Rockville, MD 20850-3173 USA

www.mesoscale.com[®]

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

Typical Standard Curve

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



Conc. (pg/mL)	sFas	
	Average Signal	%CV
0	243	6.1
1.2	401	5.1
4.9	888	4.7
20	2405	5.3
78	7719	5.3
313	29 029	8.9
1250	116 815	3.7
5000	496 984	1.7

MSD Biomarker Assays

Tested Samples

Normal human serum samples were diluted 200-fold and tested with the Human sFas Kit. Median and range of concentrations are displayed below. Concentrations are corrected for sample dilution.

Sample Type	Statistic	sFas
Serum	Median (pg/mL)	16 400
	Range (pg/mL)	10 800–2 800
	Number of Samples	20
	Samples in detectable Range	20

References

1. Papoff G, et al. An N-terminal domain shared by Fas/Apo-1 (CD95) soluble variants prevents cell death in vitro. *J. Immunol.* 1996;156:4622-4630.
2. Cascino I, et al. Soluble FAS/APO-1 Splicing Variants and Apoptosis. *Frontiers in Bioscience.* 1996 Jan 1;1:d12-18.
3. Reinehr R, et al. CD95 Ligand Is a Proliferative and Antiapoptotic Signal in Quiescent Hepatic Stellate Cells. *Gastroenterology.* 2008 May;134(5):1494-1506.
4. Del-Rey MJ, et al. Autoimmune lymphoproliferative syndrome (ALPS) in a patient with a new germline Fas gene mutation. *Immunobiology.* 2007;212(2):73-83.
5. Otsuki T, et al. Alterations of Fas and Fas-Related Molecules in Patients with Silicosis. *Exp Biol Med.* 2006 May;231(5):522-533.

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.

