

# MSD<sup>®</sup> Human TNF- $\beta$ Kit

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

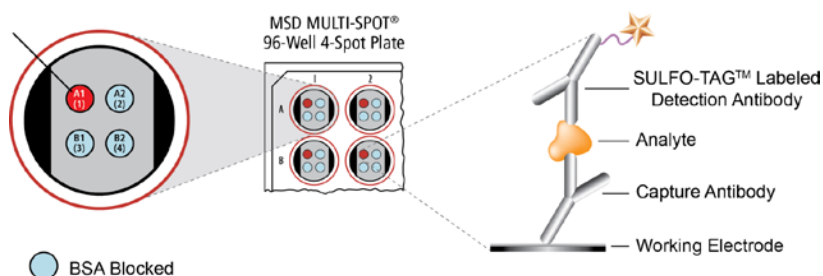


Alzheimer's Disease  
BioProcess  
Cardiac  
Cell Signaling  
Clinical Immunology  
**Cytokines**  
Growth Factors  
Hypoxia  
Immunogenicity  
Inflammation  
Metabolic  
Oncology  
Toxicology  
Vascular

## Catalog Numbers

Human TNF- $\beta$ Kit	
Kit size	
1 plate	K151LWD-1
5 plates	K151LWD-2
25 plates	K151LWD-4

## TNF- $\beta$



**Tumor necrosis factor-beta (TNF- $\beta$ )**, also known as Lymphotoxin-alpha, is a 25 kDa protein and member of the cytokine TNF-family, a group of cytokines known to contribute to apoptosis. TNF- $\beta$  is both structurally and functionally similar to TNF- $\alpha$  with 35% protein sequence homology, and both bind to the same cell surface receptors (TNF-RI and TNF-RII).<sup>1</sup> It is produced by TH1 type T-cells after antigenic or mitogenic stimulation and is cytotoxic for a range of tumor cells. Genetic polymorphism within the regulatory regions of TNF- $\alpha$  and TNF- $\beta$  (specifically 252 A to G in TNF- $\beta$ ) are linked to various cancer malignancies.<sup>2</sup> TNF- $\beta$  is a mediator of inflammatory, immunostimulatory, and antiviral responses.<sup>3</sup> It may also be involved in the proper development of secondary lymphoid organs such as lymph nodes and Peyer's patches.<sup>4</sup> TNF- $\beta$  has also been shown to contribute to the susceptibility of several autoimmune diseases, including graft-versus-host and rheumatoid arthritis<sup>5,6</sup>, as well diabetes<sup>7</sup> and other disorders.

The MSD Human TNF- $\beta$  assay is available on 96-well 4-spot plates. This datasheet outlines the performance of the assay.

## Assay Sensitivity

LLOD (pg/mL)	TNF- $\beta$
	0.23

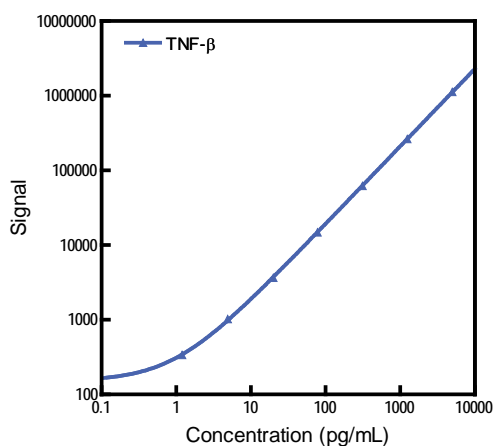
The lower limit of detection (LLOD) is a calculated concentration based on a signal 2.5 standard deviations above the background (zero calibrator blank).

## 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 TNF- $\beta$  assay.



Conc. (pg/mL)	TNF- $\beta$	
	Average Signal	%CV
0	121	6.7
1.2	339	6.0
4.9	1021	4.7
20	3642	6.6
78	14 855	6.1
313	62 095	5.1
1250	263 753	5.6
5000	1 127 795	6.0

## Company Address

MESO SCALE DISCOVERY<sup>®</sup>  
division of  
Meso Scale Diagnostics, LLC.  
1601 Research Boulevard  
Rockville, MD 20850-3173

[www.mesoscale.com](http://www.mesoscale.com)<sup>®</sup>

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

# MSD Cytokine Assays

## Specificity

The Human TNF- $\beta$  assay recognizes recombinant and native human TNF- $\beta$ . This assay shows no significant non-specific binding with following recombinant human analytes: GM-CSF, IFN- $\gamma$ , IL-1 $\beta$ , IL-2, IL-4, IL-5, IL-6, IL-8, IL-10, IL-12p40, IL-12p70, IL-13, IL-15, IL-17, TNF- $\alpha$ , or VEGF.

This assay recognizes native cynomolgus monkey TNF- $\beta$ .

## Tested Samples

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

Sample Type	Statistic	TNF- $\beta$
Serum	Median (pg/mL)	74
	Range (pg/mL)	34–96
	Number of Samples	8
	Samples in Quantitative Range	8

## References

1. Aggarwal BB, et al. Characterization of receptors for human tumour necrosis factor and their regulation by gamma-interferon. *Nature* 1985;318(6047):665-667.
2. Ibrahim A, et al. Tumor necrosis factor alpha-308 and Lymphotoxin alpha+252 genetic polymorphisms and the susceptibility to non-Hodgkin lymphoma in Egypt. *Leuk. Res.* 2012;36:694-698.
3. Remouchamps C, et al. Biology and signal transduction pathways of the Lymphotoxin- $\alpha\beta$ /LT $\beta$ R system. *Cytokine Growth Factor Rev.* 2011 Oct-Dec;22(5-6):301-10.
4. Banks TA, et al. Lymphotoxin-alpha-deficient mice: effects on secondary lymphoid organ development and humoral immune responsiveness. *J Immunol* 1995;155(4):1685-1693.
5. Chiang EY, et al. In vivo depletion of lymphotoxin-alpha expressing lymphocytes inhibits xenogeneic graft-versus-host-disease. *PLoS One.* 2012;7(3):e33106.
6. Chiang EY, et al. Targeted depletion of lymphotoxin-alpha-expressing TH1 and TH17 cells inhibits autoimmune disease. *Nat Med.* 2009 Jul;15(7):766-73.
7. Wang Y, et al. Predictive role of multilocus genetic polymorphisms in cardiovascular disease and inflammation-related genes on chronic kidney disease in Type 2 diabetes--an 8-year prospective cohort analysis of 1163 patients. *Nephrol. Dial. Transplant.* 2012;27:190-196.
8. Zhou H, et al. Collection, storage, preservation, and normalization of human urinary exosomes for biomarker discovery. *Kidney.* 2006; 69:1471-76.
9. Thomas CE, et al. Urine collection and processing for protein biomarker discovery and quantification. *Cancer Epidemiol Biomarkers & Prevention.* 2010;19:953-59.
10. Schoonenboom NS, et al. Effects of processing and storage conditions on amyloid beta (1-42) and tau concentrations in cerebrospinal fluid: implications for use in clinical practice. *Clin Chem.* 2005;51:189-95.
11. Castle JD. Purification of organelles from mammalian cells. *Curr Protoc Immunol.* 2003 Nov;Chapter 8:Unit 8.1B.

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.**

