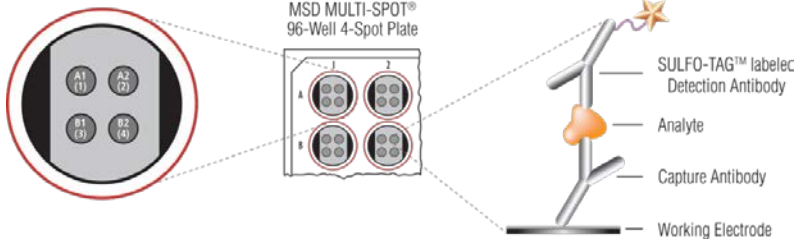


# MSD® Total IRE-1 $\alpha$ Kit

For quantitative determination in human, mouse, and rat whole cell lysate samples



1. IRE-1 $\alpha$
2. BSA blocked
3. BSA blocked
4. BSA blocked

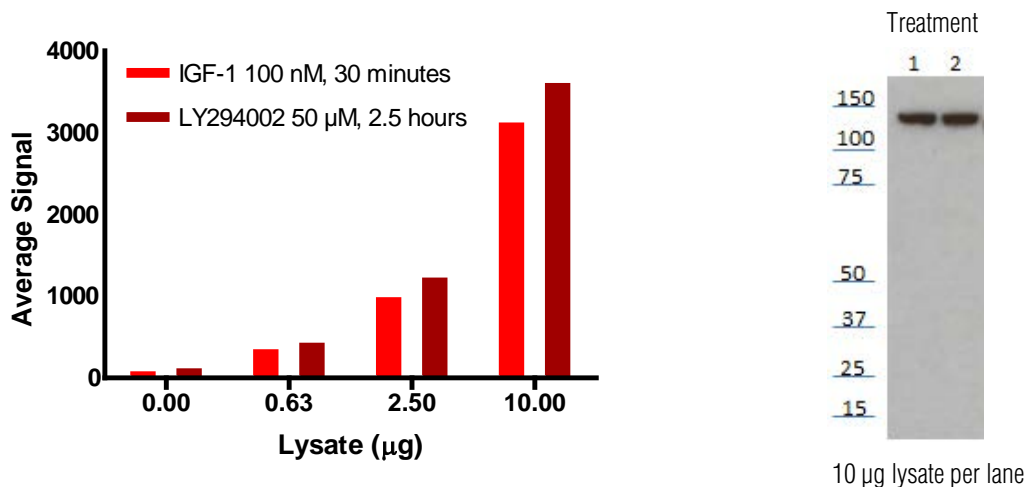


**Inositol-requiring enzyme-1 $\alpha$  (IRE-1 $\alpha$ )** is an endoplasmic reticulum (ER) transmembrane protein that mediates one of the three arms of unfolded protein response (UPR) adaptation to ER stressors. The ER is the predominant subcellular compartment for calcium storage, lipid production, and protein biosynthesis. Extracellular signaling molecules and protein receptors critical for cellular homeostasis are folded, assembled, matured, and secreted by the ER. However, ER stressors can overwhelm folding activity and unfolded proteins can accumulate in the ER, activating UPR pathways. In response to recognition of misfolded proteins in the ER lumen, IRE-1 $\alpha$  oligomerizes to activate its cytoplasmic kinase and endoribonuclease domains, which initiate cleavage of the mRNA encoding the transcription factor XBP-1 (x-box binding protein 1), an effector that drives the transcription of cytoprotective genes.<sup>1</sup> IRE-1 $\alpha$  can either return to an inactive state upon resolution of ER stress or enter a refractive state in which it no longer responds to unresolved ER stress after prolonged activation, limiting the cytoprotective duration of the UPR.<sup>2</sup> Increasing evidence suggests that ER stress and IRE-1 $\alpha$  are associated with adverse conditions such as diabetes, cancer, muscle degeneration, and neurodegenerative, bipolar, liver, cardiac, and autoimmune diseases.<sup>3-5</sup> The MSD Total IRE-1 $\alpha$  assay is available on 96-well, 4-spot plates. This datasheet outlines the performance of the assay.

## Typical Data

Representative results for the Total IRE-1 $\alpha$  Kit are illustrated below. The signal and ratio values provided are examples; individual results will vary depending upon the samples tested. Western blot analyses of each lysate type are shown for comparison.

Growing MCF-7 cells were treated with 100 nM IGF-1 for 20 minutes (treatment 1) or with 50  $\mu$ M LY294002 for 2.5 hours (treatment 2). Cell lysates were added to MSD MULTI-SPOT®, 4-spot plates coated with anti-total IRE-1 $\alpha$  antibody on one of the four spatially distinct electrodes in each well. Total IRE-1 $\alpha$  was detected with anti-total IRE-1 $\alpha$  antibody conjugated with MSD SULFO-TAG™ reagent.



**Figure 1:** Sample data generated with Total IRE-1 $\alpha$  assay. Increased signal is observed with the titration of lysates treated with 100 nM IGF-1 for 20 minutes (treatment 1) and with lysates treated with 50  $\mu$ M LY294002 for 2.5 hours (treatment 2). The Total IRE-1 $\alpha$  assay provides a quantitative measure of the data obtained with the traditional Western blot.

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

## Catalog Numbers

Total IRE-1 $\alpha$ Kit	
Kit Size	Catalog #
1 plate	K1530HD-1
5 plates	K1530HD-2
25 plates	K1530HD-4

## Ordering Information

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

## Scientific Support

Phone: 1-301-947-2025  
Email: ScientificSupport@mesoscale.com

## Company Address

MESO SCALE DISCOVERY®  
A 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 Phosphoprotein Assays

## Sample Titration

Data for positive and negative cell lysates using the Total IRE-1 $\alpha$  Kit are presented below.

Lysate ( $\mu$ g/well)	Treatment 1			Treatment 2		
	Average Signal	StdDev	%CV	Average Signal	StdDev	%CV
0	110	13	12.1	148	33	22.4
0.63	383	20	5.2	462	22	4.9
2.5	1018	18	1.7	1257	59	4.7
10	3155	100	3.2	3639	135	3.7

For a complete list of products, please visit our website at [www.mesoscale.com](http://www.mesoscale.com).

## The MSD Advantage

- **Multiplexing:** Multiple analytes can be measured in one well using typical sample volumes of 25  $\mu$ L or less without compromising speed or performance
- **Large dynamic range:** Linear range of up to five logs enables the measurement of native levels of biomarkers in normal and diseased samples without multiple dilutions
- **Minimal background:** The stimulation mechanism (electricity) is decoupled from the response (light signal), minimizing matrix interference
- **Simple protocols:** Only labels bound near the electrode surface are excited, enabling assays with fewer washes
- **Flexibility:** Labels are stable, non-radioactive, and conveniently conjugated to biological molecules

## References

1. Yoshida H, et al. XBP1 mRNA is induced by ATF6 and spliced by IRE1 in response to ER stress to produce a highly active transcription factor. *Cell*. 2001 Dec 28;107(7):881-91.
2. Li H, et al. Mammalian endoplasmic reticulum stress sensor IRE1 signals by dynamic clustering. *PNAS*. 2010 Sep 14;107(37):16113-8.
3. Wang S, et al. The impact of the unfolded protein response on human disease. *J Cell Biol*. 2012 Jun 25;197(7):857-67. Hetz C. The unfolded protein response: controlling cell fate decisions under ER stress and beyond. *Nat Rev Mol Cell Biol*. 2012 Jan 18;13(2):89-102.
4. Rath E, Haller D. Inflammation and cellular stress: a mechanistic link between immune-mediated and metabolically driven pathologies. *Eur J Nutr*. 2011 Jun;50(4):219-33.
5. Ozcan U, et al. Endoplasmic reticulum stress links obesity, insulin action, and type 2 diabetes. *Science*. 2004 Oct 15;306(5695):457-61.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR PR, SECTOR, SECTOR HTS, SULFO-TAG, [www.mesoscale.com](http://www.mesoscale.com), SMALL SPOT (design), 96 WELL 1, 4, 7, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD, MSD (design), V-PLEX, STREPTAVIDIN GOLD, and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC. © 2013 Meso Scale Diagnostics, LLC. All rights reserved.

