

MSD® CHOP Kit

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



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

Catalog Numbers

CHOP Kit	
Kit Size	Catalog #
1 plate	K150QJD-1
5 plates	K150QJD-2
25 plates	K150QJD-4

Ordering Information

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

Scientific Support

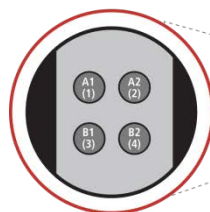
Phone: 1-301-947-2025
Email: ScientificSupport@mesoscale.com

Company Address

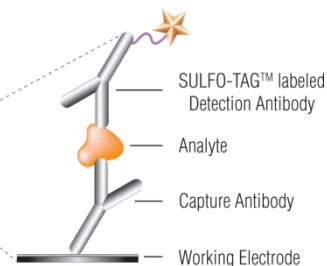
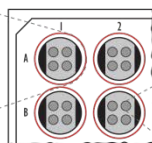
MESO SCALE DISCOVERY®
A division of
Meso Scale Diagnostics, LLC.
1601 Research Boulevard
Rockville, MD 20850-3173 USA
www.mesoscale.com®

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

1. CHOP
2. BSA blocked
3. BSA blocked
4. BSA blocked



MSD MULTI-SPOT®
96-Well 4-Spot Plate



C/EBP Homology Protein (CHOP), also known as Growth Arrest and DNA Damage Inducible Protein 153 (GADD153), DNA-damage-inducible transcript 3 (DDIT3) and C/EBP ζ , is a transcription factor that mediates one of the three arms of unfolded protein response (UPR) adaptation to endoplasmic reticulum (ER) stressors. The accumulation of unfolded or misfolded proteins in the ER is a threat to cell survival and results in a condition known as ER-stress.^{1,2} To overcome this stress, the ER initiates specific signaling pathways encompassed by the ER stress response.³ Among these are translational attenuation, upregulation of ER chaperone proteins and proteins that facilitate folding, activation of NF κ B signaling, and, as a last resort, the induction of apoptosis via CHOP.^{4,7} CHOP is a 29 kDa protein that serves as a dominant negative inhibitor of C/EBPs.⁸ CHOP is composed of an amino-terminal transcription activation domain and a carboxy-terminal basic amino-acid-rich DNA-binding domain and a leucine zipper dimerization domain. CHOP is able to act as a dominant negative inhibitor of C/EBP transcriptional activation by forming a heterodimer with C/EBPs that are normally active homodimers. The heterodimer has reduced DNA binding activity due to several key proline and glycine substitutions in the basic amino-acid-rich DNA-binding domain of CHOP. CHOP is present at low levels in the cytosol under non-stressed conditions. During ER stress, the ER stress-activated kinase PERK phosphorylates eIF2 α causing a decrease in its activity that increases the translation of ATF4 mRNA. ATF4 binds to and activates the CHOP promoter thereby increasing its levels during ER stress.⁹ While the precise role of CHOP in ER stress-induced apoptosis is not completely understood, one of the most widely cited mechanisms is via the suppression of the pro-survival protein Bcl-2.¹⁰ CHOP is involved in the response of many diseases (ischemia, neurodegeneration, cancer, diabetes) which makes this an attractive marker for many therapeutic areas. The MSD CHOP assay is available on 96-well, 4-spot plates. This datasheet outlines the performance of the assay.

Typical Data

Representative results for the CHOP 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.

This assay was developed using recombinant human CHOP protein (data shown below) as well as with the nuclear fraction of rat hepatoma cells treated with ER stress-inducing agents, thapsigargin and tunicamycin (data not shown). Recombinant protein or nuclear cell lysate fractions were added to MSD MULTI-SPOT 4-spot plates coated with anti-CHOP antibody on one of the four spatially distinct electrodes in each well. CHOP was detected with anti-CHOP antibody conjugated with MSD SULFO-TAG reagent.

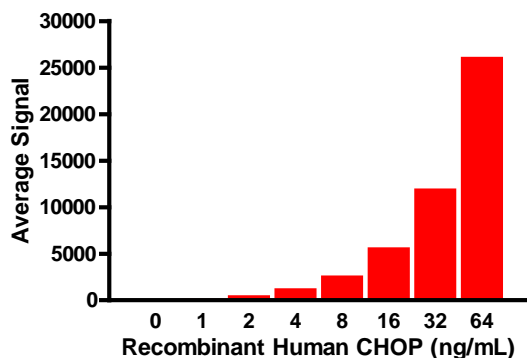


Figure 1: Sample data generated with CHOP assay. Increased signal is observed with the titration of recombinant human CHOP protein.

MSD Phosphoprotein Assays

Sample Titration

Data for recombinant CHOP protein using the CHOP Kit are presented below.

Sample (ng/well)	CHOP recombinant protein		
	Average Signal	StdDev	%CV
0	52	2	4.1
1.0	399	13	3.2
2.0	754	21	2.8
4.0	1509	37	2.5
8.0	2888	150	5.2
16	5900	122	2.1
32	12244	332	2.7
64	26383	268	1.0

For a complete list of products, please visit our website at www.mesoscale.com.

The MSD Advantage

- **Multiplexing:** Multiple analytes can be measured in one well using typical sample volumes of 25 μ 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. Kopito RR. Aggresomes, inclusion bodies and protein aggregation. *Trends Cell Biol.* 2000 Dec;10(12):524-30.
2. Kaufman RJ. Stress signaling from the lumen of the endoplasmic reticulum: coordination of gene transcriptional and translational controls. *Genes Dev.* 1999 May 15;13(10):1211-33.
3. Mori K. Tripartite management of unfolded proteins in the endoplasmic reticulum *Cell.* 2000 May 26;101(5):451-4.
4. Harding HP, et al. Transcriptional and translational control in the Mammalian unfolded protein response. *Annu Rev Cell Dev Biol.* 2002;18:575-99.
5. Kozutsumi MY, et al. The presence of malfolded proteins in the endoplasmic reticulum signals the induction of glucose-regulated proteins. *Nature.* 1988 Mar 31;332(6163):462-4.
6. Pahl HL, et al. Activation of transcription factor NF-kappaB by the adenovirus E3/19K protein requires its ER retention. *J Cell Biol.* 1996 Feb;132(4):511-22.
7. Matsumoto M, et al. Ectopic expression of CHOP (GADD153) induces apoptosis in M1 myeloblastic leukemia cells. *FEBS Lett.* 1996 Oct 21;395(2-3):143-7.
8. Ron D, Habener JF. CHOP, a novel developmentally regulated nuclear protein that dimerizes with transcription factors C/EBP and LAP and functions as a dominant-negative inhibitor of gene transcription. *Genes Dev.* 1992 Mar;6(3):439-53.
9. Wang XZ, et al. Signals from the stressed endoplasmic reticulum induce C/EBP-homologous protein (CHOP/GADD153). *Mol Cell Biol.* 1996 Aug;16(8):4273-80.
10. McCullough KD, et al. Gadd153 sensitizes cells to endoplasmic reticulum stress by down-regulating Bcl2 and perturbing the cellular redox state. *Mol Cell Biol.* 2001 Feb;21(4):1249-59.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR PR, SECTOR, SECTOR HTS, SULFO-TAG, 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.

