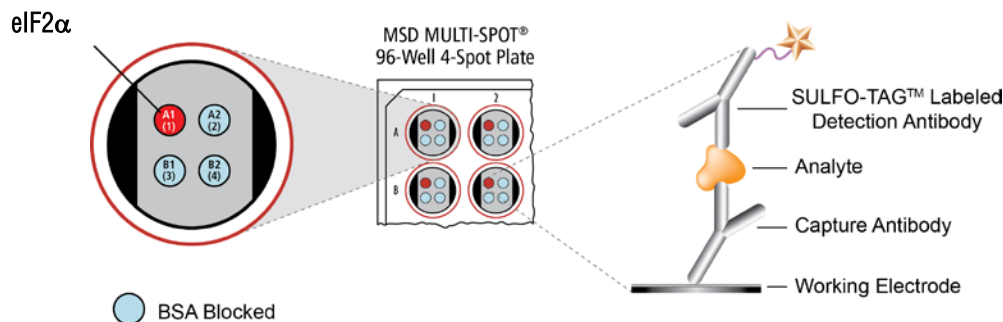


MSD® Total eIF2 α Whole Cell Lysate Kit

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



Eukaryotic protein synthesis is a tightly coordinated process consisting of initiation, elongation, and termination phases. Initiation of mRNA translation requires binding of Met-tRNA to the 40S ribosomal subunit.

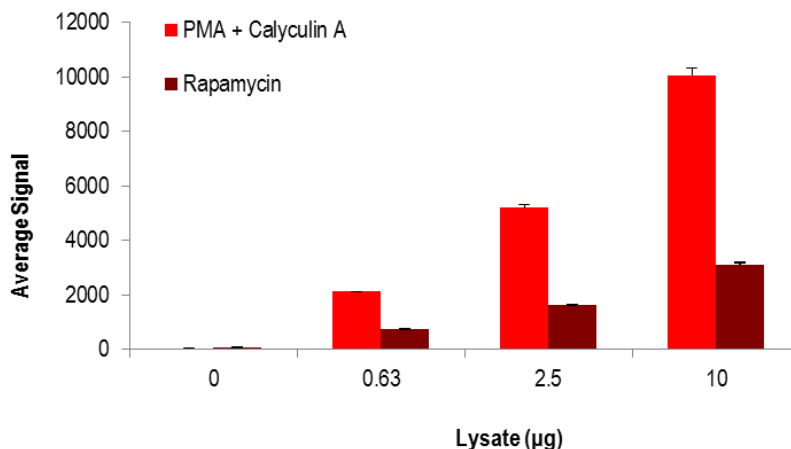
Eukaryotic translation initiation factor 2 (eIF2), a heterodimer consisting of alpha, beta, and gamma subunits, is an essential regulator of translational initiation. Active eIF2-GTP, Met-tRNA, and the 40S subunit form the ternary 43S preinitiation complex.^{1,2} Once initiation is completed, inactive eIF2-GDP complex is ejected from the ribosome, and the guanine nucleotide exchange reaction catalyzed by eIF2 β recycles eIF2 to an active state, permitting additional rounds of initiation.^{1,2} The alpha subunit, eIF2 α , is the regulatory domain of the eIF2 heterotrimer; phosphorylation of eIF2 α (Ser51) stabilizes the eIF2-GDP-eIF2 β complex, globally repressing translation.^{1,2}

Four eIF2 α kinases have been identified to date, and each kinase responds to a different stressor. General control non-depressible-2 (GCN2) is activated during amino acid starvation;³ protein kinase R (PKR) is activated in response to dsRNA;⁴ PKR-like endoplasmic reticulum kinase is activated in response to accumulation of misfolded proteins in the endoplasmic reticulum;⁵ and heme-regulated inhibitor (HRI) limits protein synthesis in heme-deficient cells.⁶

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

Typical Data

Representative results for the kit are illustrated below. The signal and ratio values provided are examples; individual results may vary depending upon the samples tested. with MSD SULFO-TAG reagent.



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

Catalog Numbers

Total eIF2 α
Whole Cell Lysate Kit

Kit Size

1 plate K150NGD-1

5 plates K150NGD-2

25 plates K150NGD-4

Ordering information

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

Lysate Titration

Data for each treatment using the [Meso Scale Discovery PhosphoScan[®]](#) Kit is presented below.

Lysate ($\mu\text{g}/\text{Well}$)	200 nM PMA+50 nM Calyculin A			1 μM Rapamycin		
	Average Signal	StdDev	%CV	Average Signal	StdDev	%CV
0	42	5	11.9	57	6	9.9
0.63	2119	19	0.9	736	5	0.7
2.5	5196	122	2.3	1633	7	0.4
10	10060	273	2.7	3118	88	2.8

MSD Advantage

- **Multiplexing:** Multiple analytes can be measured in one well using typical sample amounts of 25 $\mu\text{g}/\text{well}$ 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 signal (light)
- **Simple protocols:** Only labels near the electrode surface are detected, enabling no-wash assays
- **Flexibility:** Labels are stable, non-radioactive, and conveniently conjugated to biological molecules
- **High sensitivity and precision:** Multiple excitation cycles of each label enhance light levels and improve sensitivity

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

References

1. Hershey JW. Protein phosphorylation controls translation rates. *J Biol Chem.* 1989 Dec 15; 264(35):20823-6.
2. Kimball SR. Eukaryotic initiation factor eIF2. *Int J Biochem Cell Biol.* 1999 Jan;31(1):25-29.
3. Grallert B, Boye E. The Gcn2 kinase as a cell cycle regulator. *Cell Cycle.* 2007 Nov 15;6(22):2768-72.
4. Nalagatla SR, et al. Regulation of innate immunity through RNA structure and the protein kinase PKR. *Curr Opin Struct Biol.* 2011 Feb; 21(1):119-27.
5. Raven JF, Koromilas AE. PERK and PKR: Old kinases learn new tricks. *Cell Cycle.* 2008 May 1;7(9):1146-50.
6. Chen JJ. Regulation of protein synthesis by the heme regulated eIF2alpha kinase: relevance to anemias. *Blood.* 2007 Apr 1;109(7):2693-9.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, WWW.MESOSCALE.COM, MSD, MSD (DESIGN), DISCOVERY WORKBENCH, QUICKPLEX, MULTI-ARRAY, MULTI-SPOT, SULFO-TAG, 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.

