

MSD® Phospho(Ser112)/Total BAD Assay Whole Cell Lysate Kit

For quantitative determination in human and monkey whole cell lysate samples



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

Catalog Numbers

Phospho(Ser112)/Total BAD Assay: Whole Cell Lysate Kit

Kit size

1 plate	K15103D-1
5 plates	K15103D-2
20 plates	K15103D-3

Phospho-BAD Whole Cell Lysate Set	
200 µg	C11CC-1

Ordering information

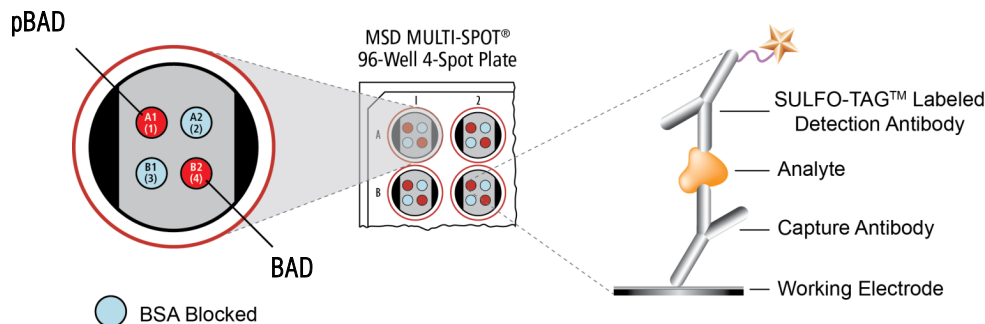
MSD Customer Service
Phone: 1-301-947-2085
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com

Company Address

MESO SCALE DISCOVERY®
A division of
Meso Scale Diagnostics, LLC.
9238 Gaither Road
Gaithersburg, MD 20877 USA

www.mesoscale.com®

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



Bcl-2-antagonist of cell death protein (BAD), a member of the pro-apoptotic Bcl-2 family of proteins, functions by displacing the binding of Bax to Bcl-2 and Bcl-xL, and causes cell death by apoptosis. The binding of cytokines and growth factors to cell surface receptors activates intracellular signal transduction cascades that promote cell survival. Akt phosphorylates BAD on Ser136. BAD is also phosphorylated by protein kinase A (PKA) and p90 ribosomal S6 kinase (p90RSK) on Ser112. Phosphorylated BAD binds to members of the 14-3-3 protein family. This inhibits its interaction with Bcl-2 and Bcl-xL through cytosolic sequestration. The anti-apoptotic proteins Bcl-2 and Bcl-xL are then free to interact with Apaf-1 and BID, thus promoting cell survival.

The MSD Phospho(Ser112)/Total BAD Assay is available on 96-well 4-Spot plates. This datasheet outlines the performance of the assay.

Typical Data

Representative results for the Phospho(Ser112)/Total BAD Assay are illustrated below. The signal and ratio values provided below are example data; individual results may vary depending upon the samples tested. Western blot analyses of each lysate type were performed with phospho-BAD (Ser112) and total BAD antibodies and are shown below for comparison.

Serum deprived COS-7 cells were treated with staurosporine (1 mM, 3 hours) (negative), or treated with PMA (200 nM, 1 hour) (positive). Whole cell lysates were added to MSD MULTI-SPOT® 4-Spot plates coated with anti-phospho-BAD (Ser112) antibody and anti-total BAD antibody on spatially distinct electrodes within a well. Phosphorylated and total BAD were detected with anti-total BAD antibody conjugated with MSD SULFO-TAG™ reagent.

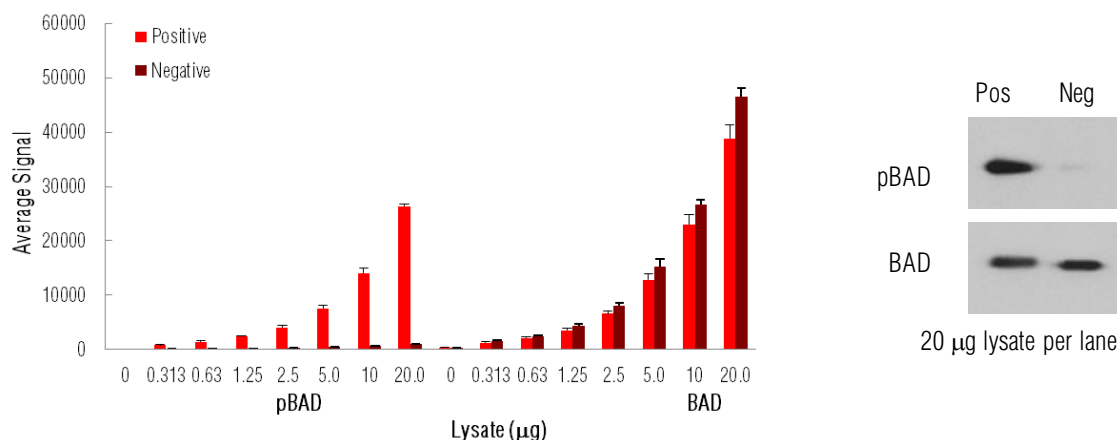


Fig. 1: Sample data generated with the MULTI-SPOT Phospho(Ser112)/Total BAD Assay. Increased signal for phosphorylated BAD was observed with only pBAD positive cell lysate. Total BAD signal increased throughout the titration of both pBAD positive and negative cell lysates. The Phospho(Ser112)/Total BAD Assay provides a quantitative measure of the data obtained with the traditional Western blot.

MSD Phosphoprotein Assays

Lysate Titration

Data for pBAD positive and negative COS-7 cell lysates using the MULTI-SPOT Phospho(Ser112)/Total BAD Assay are presented below.

	Lysate (μ g)	Positive			Negative			P/N
		Average Signal	StdDev	%CV	Average Signal	StdDev	%CV	
pBAD	0	22	5	23.0	21	8	37.9	
	0.313	807	82	10.2	92	4	3.9	8.8
	0.63	1396	151	10.8	133	6	4.8	11
	1.25	2466	110	4.5	179	7	4.0	14
	2.5	4051	301	7.4	248	4	1.4	16
	5.0	7547	599	7.9	426	5	1.2	18
	10	14099	906	6.4	666	2	0.3	21
	20.0	26366	395	1.5	988	5	0.5	27
BAD	0	321	31	9.7	349	18	5.1	
	0.313	1260	139	11.0	1488	144	9.7	0.8
	0.63	2077	206	9.9	2435	206	8.5	0.9
	1.25	3532	427	12.1	4360	303	6.9	0.8
	2.5	6552	556	8.5	8050	519	6.4	0.8
	5.0	12852	1125	8.8	15246	1343	8.8	0.8
	10	23059	1794	7.8	26751	789	2.9	0.9
	20.0	38839	2518	6.5	46542	1486	3.2	0.8

MSD Advantage

- **Multiplexing:** Multiple analytes can be measured in one well using typical sample amounts of 25 μ g/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

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.
© 2011 Meso Scale Diagnostics, LLC. All rights reserved.

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

