

MSD[®] MULTI-ARRAY Assay System

Human DJ-1/PARK7 Kit

1-Plate Kit	K151THD-1
5-Plate Kit	K151THD-2
25-Plate Kit	K151THD-4



MSD Neurodegenerative Disease Assays

Human DJ-1/PARK7 Kits

For use with human cerebrospinal fluid (CSF) and serum.

This package insert must be read in its entirety before using this product.

FOR RESEARCH USE ONLY.

NOT FOR USE IN DIAGNOSTIC PROCEDURES.

MESO SCALE DISCOVERY[®]

A division of Meso Scale Diagnostics, LLC.

1601 Research Boulevard

Rockville, MD 20850-3173 USA

www.mesoscale.com

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

Table of Contents

Introduction	4
Principle of the Assay	4
Additional Materials and Equipment.....	5
Safety	6
Best Practices and Technical Hints	6
Reagent Preparation	7
Protocol	9
Curve Fitting.....	9
Typical Data.....	10
Sensitivity	10
Assay Components	10
References.....	11
Summary Protocol	13
Plate Diagram	15

Ordering Information

MSD Customer Service

Phone: 1-301-947-2085, 1-240-632-5885
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com

MSD Scientific Support

Phone: 1-301-947-2025
Fax: 1-240-632-2219 attn: Scientific Support
Email: ScientificSupport@mesoscale.com

Introduction

The DJ-1/PARK7 gene is linked to familial early-onset Parkinson's disease (PD) and encodes a 189 amino acid protein belonging to the DJ-1/Thi/Pfpl superfamily of proteins.¹ DJ-1/PARK7 has been implicated as an oxidative stress sensor that functions to promote the survival of dopaminergic neurons.¹ Under oxidative conditions, DJ-1/PARK7 undergoes a pI shift from 6.2 to 5.8 and protects against mitochondrial damage.²⁻⁴ DJ-1/PARK7 also functions as a redox-dependent protein chaperone to suppress aggregation of α -synuclein, the main constituent in Lewy bodies prominent in the brains of PD patients.⁵ Furthermore, DJ-1/PARK7 forms a complex with parkin and PTEN-induced putative kinase 1 (PINK1) during oxidative stress to promote the degradation of parkin substrates, many of which have been implicated in neurodegeneration.^{6,7} Along with α -synuclein, DJ-1 has emerged as a candidate biomarker of PD in human CSF. The MSD Human DJ-1/PARK7 Kit can be used to measure DJ-1/PARK7 in human CSF and serum.

Principle of the Assay

MSD neurodegenerative disease assays provide a rapid and convenient method for measuring the levels of protein targets within a single, small-volume sample. Human DJ-1/PARK7 is a sandwich immunoassay. MSD provides a plate pre-coated with capture antibodies on independent and well-defined spots in the layout shown below. The user adds the sample and a solution containing detection antibodies conjugated with electrochemiluminescent labels (MSD SULFO-TAG™) over the course of one or more incubation periods. Analytes in the sample bind to capture antibodies immobilized on the working electrode surface; recruitment of the detection antibodies by the bound analytes completes the sandwich. The user adds an MSD buffer that provides the appropriate chemical environment for electrochemiluminescence and loads the plate into an MSD instrument where a voltage applied to the plate electrodes causes the captured labels to emit light. The instrument measures the intensity of emitted light to provide a quantitative measure of analyte in the sample.

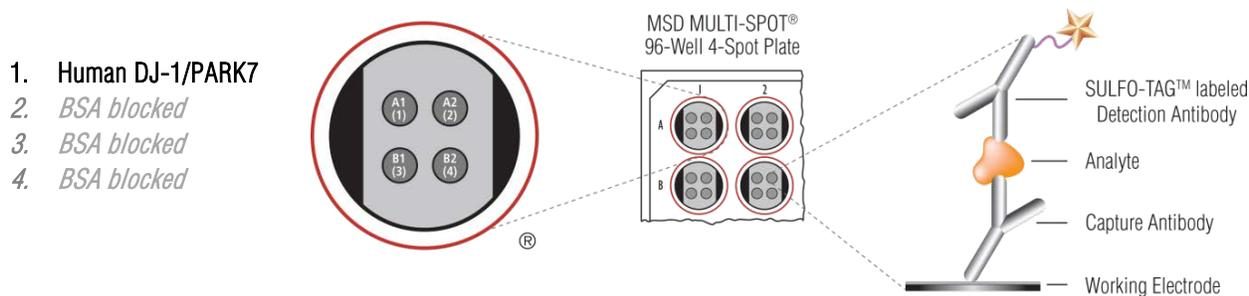


Figure 1. Spot diagram showing placement of analyte capture antibodies for the Human DJ-1/PARK7 Kit. The numbering convention for the different spots is maintained in the software visualization tools, on the plate packaging, and in the data files.

Reagents Supplied

Reagent	Storage	Catalog #	Size	Quantity Supplied			Description
				1 Plate Kit	5 Plate Kit	25 Plate Kit	
MULTI-SPOT® 96-well 4-spot Human DJ-1 Plate	2–8°C	N451THA-1	4-Spot	1	5	25	96-well plate, foil sealed with desiccant.
SULFO-TAG Anti- Human DJ-1 Antibody (50X) ¹	2–8°C	D21TH-2	75 µL	1 vial			SULFO-TAG–conjugated antibody
		D21TH-3	375 µL		1 vial	5 vials	
Human DJ-1 Calibrator (20X) ²	≤-70°C	C01TH-2	30 µL	1 vial	5 vials	25 vials	Purified recombinant human DJ-/PARK7 protein in a buffered protein diluent.
Diluent 35	2–8°C	R50AE-3	30 mL	1 bottle			Diluent for samples and calibrator; contains blockers and preservatives.
		R50AE-2	150 mL		1 bottle	5 bottles	
Diluent 100	2–8°C	R50AA-4	50 mL	1 bottle	1 bottle	5 bottles	Diluent for detection antibody; contains protein, blockers, and preservatives.
Blocker D–G ³	≤-10°C	R93BH-3	1.0 mL	1 vial	2 vials	10 vials	Goat gamma globulin solution
Blocker D–M ³	≤-10°C	R93BM-1	200 µL	1 vial			Mouse gamma-globulin solution
		R93BM-2	900 µL		1 vial	5 vials	
Blocker D–R ³	≤-10°C	R93BR-1	50 µL	1 vial			Rabbit gamma-globulin solution
		R93BR-2	200 µL		1 vial	5 vials	
Read Buffer T (4X)	RT	R92TC-3	50 mL	1 bottle	1 bottle	5 bottles	MSD buffer to catalyze the electro-chemiluminescence reaction

Additional Materials and Equipment

- Appropriately sized tubes for reagent preparation
- Polypropylene microcentrifuge tubes for preparing dilutions
- Liquid handling equipment for desired throughput, capable of dispensing 10 to 150 µL/well into a 96-well microtiter plate
- Plate washing equipment: automated plate washer or multichannel pipette
- Microtiter plate shaker (rotary) capable of shaking at 300–1000 rpm.
- Phosphate-buffered saline plus 0.05% Tween-20 for plate washing or MSD Wash Buffer, catalog # R61AA-1
- Adhesive plate seals
- Deionized water

¹ SULFO-TAG conjugated detection antibodies should be stored in the dark.

² **Biohazard Statement:** This product was derived from human material and should be considered potentially infectious. Appropriate precautions should be used when handling this material.

³ Blockers D–G, D–M, and D–R can tolerate up to 5 freeze–thaw cycles. Alternatively, aliquots of Blockers D–G, D–M, and D–R can be stored at 2–8°C for up to 1 month.

Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling kit components. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines.

Additional product-specific safety information is available in the safety data sheet, which can be obtained from MSD Customer Service.

Best Practices and Technical Hints

- Do not mix or substitute reagents from different sources or different kit lots.
- Dilute calibrators and samples in polypropylene microcentrifuge tubes; use a fresh pipette tip for each dilution; vortex after each dilution before proceeding.
- Avoid prolonged exposure of detection antibody (stock or diluted) to light. During the antibody incubation step, plates do not need to be shielded from light except for direct sunlight.
- Shaking should be vigorous with a rotary motion between 300 and 1000 rpm.
- Avoid bubbles in wells at all pipetting steps. Bubbles may lead to variable results; bubbles introduced when adding read buffer may interfere with signal detection. Do not shake plate after adding read buffer.
- Use reverse pipetting when necessary to avoid introduction of bubbles, and pipette to the bottom corner of empty wells.
- When using an automated plate washer, rotating the plate 180 degrees between wash steps may improve assay precision.
- Gently tap the plate to remove residual fluid after washing.
- Read buffer should be at room temperature when added to the plate.
- Keeping time intervals consistent between adding read buffer and reading the plate should improve inter-plate precision. Limit the time the plate is incubated with read buffer.
- Remove plate seals prior to reading the plate.
- If an incubation step needs to be extended, avoid letting the plate dry out by keeping sample or detection antibody solution in the wells.
- If assay results are above the top of the calibration curve, dilute samples, and repeat the assay.
- When running partial plates, use the sector map in the instrument or software manual to select the wells to be used. Seal the unused portion of the plate with a plate seal to avoid contaminating unused wells. After reading a partial plate, remove fluid, reseal unused sectors, return plate to its original foil pouch with desiccant pack, and seal pouch with tape. Partially used plates may be stored for up to 14 days at 2–8°C.
- You may adjust volumes proportionally when preparing reagents.

Reagent Preparation

Bring all reagents to room temperature.

Prepare Calibrator Dilutions

MSD supplies calibrator for the Human DJ-1/PARK7 Kit at a concentration that is 20-fold higher than the recommended highest standard. We recommend a 7-point calibration curve with 4-fold serial dilution steps and a zero calibrator blank. Thaw the stock calibrator and keep on ice, then add to diluent at room temperature to make the calibration curve solutions.

Calibrator	Human DJ-1/PARK7 (pg/mL)	Dilution Factor
Stock Calibrator	2 000 000	
Calibrator-01	100 000	20
Calibrator-02	25 000	4
Calibrator-03	6 250	4
Calibrator-04	1 563	4
Calibrator-05	391	4
Calibrator-06	98	4
Calibrator-07	24	4
Calibrator-08	0	n/a

To prepare 7 calibration solutions plus a zero calibrator blank for up to 4 replicates:

1. Prepare the highest calibrator by adding 15 μ L of stock calibrator to 285 μ L of Diluent 35. Mix well.
2. Prepare the next calibrator by transferring 50 μ L of the highest calibrator to 150 μ L of Diluent 35. Mix well. Repeat 4-fold serial dilution 5 times to generate 7 calibrators.
3. Use Diluent 35 as the blank.

Sample Collection and Handling

CSF

Sample collection methods and pre-analytical conditions may cause variability in measured analyte levels.⁸⁻¹⁰ MSD recommends reviewing current literature and protocols for collection and handling of CSF samples.

Serum and Plasma

Plasma prepared in heparin tubes commonly displays additional clotting following thawing of the sample. Remove clots and all solid material by centrifugation. Avoid multiple freeze–thaw cycles for serum and plasma samples.

Dilute Samples

Dilute samples with Diluent 35. For human CSF and serum samples, MSD recommends a minimum 8-fold sample dilution; however depending on the sample set under investigation, you may need to use a higher dilution factor. For example, to dilute 8-fold, add 20 μ L of sample to 140 μ L of Diluent 35.

Prepare Detection Antibody Solution

MSD provides detection antibody as a 50X stock solution. The working solution is 1X. Prepare the detection antibody solution immediately prior to use.

For 1 plate, combine:

- 60 μ L of 50X SULFO-TAG Anti-Human DJ-1/PARK7 Antibody
- 300 μ L of 10% Blocker D-G (1% final concentration in detection antibody solution)
- 150 μ L of 2% Blocker D-M (0.1% final concentration in detection antibody solution)
- 30 μ L of 10% Blocker D-R (0.1% final concentration in detection antibody solution)
- 2460 μ L of Diluent 100

Prepare Read Buffer

MSD provides Read Buffer T as a 4X stock solution. The working solution is 2X.

For 1 plate, combine:

- 10 mL of Read Buffer T (4X)
- 10 mL of deionized water

You may keep excess diluted read buffer in a tightly sealed container at room temperature for up to 1 month.

Prepare MSD Plate

MSD plates are pre-coated with capture antibodies (Figure 1) and exposed to a proprietary stabilizing treatment to ensure the integrity and stability of the immobilized antibodies. Plates may be used as delivered; no additional preparation (e.g., pre-wetting) is required.

Protocol

1. **Block plate.** Add 150 μL of Diluent 35 to each well. Seal the plate with an adhesive plate seal and incubate at room temperature with shaking for 1 hour.

You may prepare calibrators, samples, and detection antibody during incubation.

2. **Wash, Add Detection Antibody Solution and Sample:** Wash the plate 3 times with at least 150 μL /well of PBS-T. Add 25 μL of detection antibody solution to each well. Add 25 μL of diluted sample or calibrator per well. Seal the plate with an adhesive plate seal and incubate at room temperature with shaking for 2 hours.

You may prepare diluted read buffer during incubation.

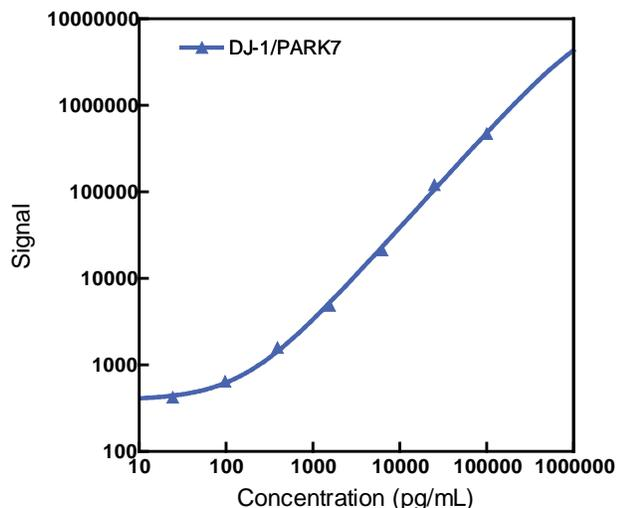
3. **Wash and Read:** Wash the plate 3 times with at least 150 μL /well of PBS-T. Add 150 μL of 2X Read Buffer T to each well. Read plate on the MSD instrument. No incubation in read buffer is required before reading the plate.

Curve Fitting

Run at least 1 set of calibrators in duplicate to generate the calibration curve. The calibration curve is modeled using least squares fitting algorithms so that signals from the calibrators can be used to calculate the concentration of analyte in the samples. The assay has a wide dynamic range (4 logs), which allows for accurate quantification in samples without the need for multiple dilutions or repeated testing. The data displayed below were generated by DISCOVERY WORKBENCH[®] analysis software using a 4-parameter, logistic curve-fitting model (sigmoidal dose-response) with a $1/Y^2$ weighting function, which provides a better fit of data over a wide dynamic range, particularly at the low end of the calibration curve.

Typical Data

The following calibration curve graph illustrates the dynamic range of the assay. Actual signals will vary. Best quantification of unknown samples will be achieved by generating a calibration curve for each plate using a minimum of 2 replicates of calibrators.



Human DJ-1/PARK7		
Conc. (pg/mL)	Average Signal	%CV
0	321	0.9
24	422	4.4
98	646	1.6
391	1 595	1.8
1 563	4 874	0.6
6 250	21 319	0.8
25 000	120 468	1.2
100 000	471 893	2.2

Sensitivity

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

Human DJ-1/PARK7	
Average LLOD (pg/mL)	12.0

Assay Components

Calibrators

The assay calibrator uses purified human DJ-1/PARK7.

Antibodies

Analyte	Source Species		Assay Generation
	MSD Capture Antibody	MSD Detection Antibody	
Human DJ-1/PARK7	Rat Monoclonal	Goat Polyclonal	A

References

1. Thomas B, et al. Parkinson's disease. *Human Molecular Genetics*. 2007;16:R183-4.
2. Canet-Aviles RM, et al. The Parkinson's disease protein DJ-1 is neuroprotective due to cysteine-sulfinic acid-driven mitochondrial localization. *PNAS*. 2004;101:9103-8.
3. Mitsumoto A, et al. Oxidized forms of peroxiredoxins and DJ-1 on two-dimensional gels increased in response to sublethal levels of paraquat. *Free Radical Res*. 2001;35:301-10.
4. Mitsumoto A, et al. DJ-1 is an indicator for endogenous reactive oxygen species elicited by endotoxin. *Free Radical Res*. 2001;35:885-93.
5. Shendelman S, et al. DJ-1 is a redox-dependent molecular chaperone that inhibits α -synuclein aggregate formation. *PLoS Biol*. 2004;2:1764-73.
6. Dawson TM. Parkin and defective ubiquitination in Parkinson's disease. *J. Neural Transm. Suppl*. 2006;70:209-13.
7. Moore DJ, et al. Association of DJ-1 and parkin mediated by pathogenic DJ-1 mutations and oxidative stress. *Hum Mol Genet* 2005;14:71-84.
8. Shi M, et al. Significance and confounders of peripheral DJ-1 and alpha-synuclein in Parkinson disease. *Neurosci Lett*. 2010;480(1):78-82.
9. Mattsson N, et al. Inter-laboratory variation in cerebrospinal fluid biomarkers for Alzheimer's disease: united we stand, divided we fall. *Clin Chem Lab Med*. 2010;48:603-7.
10. Zetterberg H, et al. Clinical proteomics in neurodegenerative disorders. *Acta Neurol Scand*. 2008;118:1-11.

Summary Protocol

Human DJ-1/PARK7 Kits

*MSD provides this summary protocol for your convenience.
Please read the entire detailed protocol prior to performing
the Human DJ-1/PARK7 assays.*

Sample and Reagent Preparation

Bring all reagents to room temperature. The calibration curve, diluted samples, and detection antibody solution should be prepared during step 1 and used within one hour of preparation.

Prepare 7 calibration solutions in Diluent 35 using the supplied calibrator:

- Dilute the stock calibrator by adding 15 μL of stock calibrator to 285 μL of Diluent 35. Mix well.
- Perform a series of 4-fold dilution steps and prepare a zero calibrator blank.

Dilute samples 8-fold in Diluent 35 before adding to the plate.

Prepare detection antibody solution by diluting stock detection antibody and blockers in Diluent 100.

Prepare 2X Read Buffer T by diluting stock 4X Read Buffer T 2-fold with deionized water.

Step 1: Block Plate

Add 150 μL /well of Diluent 35.

Incubate at room temperature with shaking for 1 hour.

Step 2: Wash, Add Detection Antibody Solution and Sample

Wash plate 3 times with at least 150 μL /well of PBS-T.

Add 25 μL /well of 1X detection antibody solution.

Add 25 μL /well of sample (calibrators or diluted samples).

Incubate at room temperature with shaking for 2 hours.

Step 3: Wash and Read Plate

Wash plate 3 times with at least 150 μL /well of PBS-T.

Add 150 μL /well of 2X Read Buffer T.

Analyze plate on MSD instrument.

Plate Diagram

