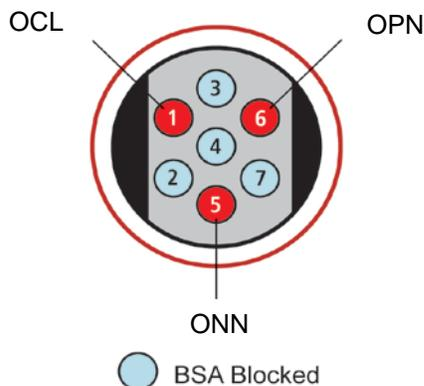


# MSD<sup>®</sup> 96-Well MULTI-SPOT<sup>®</sup> Human Bone Panel II

The following assay protocol has been optimized for analysis of osteocalcin (OCL), osteonectin (ONN), and osteopontin (OPN) in human serum and plasma samples.

	Storage
<b>MSD Materials</b>	
<input type="checkbox"/> MULTI-SPOT 96-well 7-Spot Human Bone Panel II Plate(s)	2–8°C
<input type="checkbox"/> SULFO-TAG <sup>™</sup> Anti-hOsteocalcin Antibody (50X) <sup>1</sup>	2–8°C
<input type="checkbox"/> SULFO-TAG Anti-hOsteonectin Antibody (50X) <sup>1</sup>	2–8°C
<input type="checkbox"/> SULFO-TAG Anti-hOsteopontin Antibody (50X) <sup>1</sup>	2–8°C
<input type="checkbox"/> Human Bone Panel II Calibrator Blend Osteocalcin (OCL – 0.2 µg/mL) Osteonectin <sup>2</sup> (ONN – 2 µg/mL) Osteopontin (OPN – 0.2 µg/mL)	≤-70°C
<input type="checkbox"/> Diluent 7	≤-10°C
<input type="checkbox"/> Diluent 11	≤-10°C
<input type="checkbox"/> Blocker A Kit	RT
<input type="checkbox"/> MSD Read Buffer T (4X)	RT



The SECTOR<sup>®</sup> Imager data file will identify spots according to their well location, not by the coated capture antibody name.

<sup>1</sup> SULFO-TAG-conjugated detection antibodies should be stored in the dark.

<sup>2</sup> Osteonectin in the Calibrator Blend is derived from human source material which has been tested and found to be negative or non reactive for HBsAg, anti-HB core, Syphilis, anti-HVC, anti- HIV-1 antigens, anti-HIV1/2 and anti-HTLV1/2. This material should be handled and disposed of in accordance with local, state, and federal guidelines.

FOR RESEARCH USE ONLY.  
NOT FOR USE IN DIAGNOSTIC PROCEDURES.



**Note:** 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 (SDS), which can be obtained from MSD Customer Service.

**Notes:**

## *Other Materials & Equipment (not supplied)*

- Deionized water for diluting concentrated buffers
- Phosphate buffered saline (PBS) for plate washing
- 15 mL and 50 mL tubes
- Adhesive plate seals
- Microtiter plate shaker
- Microcentrifuge tubes for making dilutions of standards
- Automated plate washer or other efficient multi-channel pipetting equipment for washing 96-well plates
- Appropriate liquid handling equipment for desired throughput capable of accurately dispensing 25  $\mu$ L and 150  $\mu$ L into a 96-well microplate

## *Protocol at a Glance*

- Step 1.** Add Blocking Solution, incubate 1 hour, wash.
- Step 2.** Add 25  $\mu$ L of Diluent 7.  
Add 25  $\mu$ L of Calibrator or diluted samples (diluted 20-fold), incubate 2 hours, wash.
- Step 3.** Add 25  $\mu$ L of Detection Antibody, incubate 1 hour, wash.
- Step 4.** Add 150  $\mu$ L of Read Buffer, read plate, and analyze data.

*Read the entire detailed instructions before beginning work.*

The full protocol that follows describes the most conservative approach to achieving highly sensitive results using MSD technology to quantify OCL, ONN, and OPN. The protocol can be completed in approximately 4.5 hours if each reagent is prepared during the preceding incubation. This time can be reduced to **3.5 hours if the blocking step is performed overnight** prior to performing the assay.



## Preparation Instructions

**Thaw Reagents.** Thaw MSD Human Bone Panel II Calibrator Blend, Diluent 7, and Diluent 11. Vortex briefly once thawed. If there is a precipitate, mix gently, and warm to room temperature to dissolve. Keep all materials on ice until use. The remaining amounts of the diluents after use can be aliquoted and refrozen as needed.

**Prepare Blocker A Solution.** Prepare MSD Blocker A solution following the instructions included in the MSD Blocker A Kit.

**Prepare Detection Antibody Solution.** MSD provides the detection antibodies blended in a 50X stock solution. The working detection antibody solution is 1X. You will need 3 mL per plate.

In a 15 mL tube, combine:

- 2.82 mL Diluent 11
- 60 µL of 50X SULFO-TAG Anti-hOsteocalcin Antibody
- 60 µL of 50X SULFO-TAG Anti-hOsteonectin Antibody
- 60 µL of 50X SULFO-TAG Anti-hOsteopontin Antibody

**Prepare Read Buffer T.** MSD provides Read Buffer T as a 4X stock solution. The working solution is 1X. You will need 20 mL per plate at a final 1X concentration.

In a 50 mL tube, combine:

- 15 mL deionized water
- 5 mL 4X Read Buffer T

### Prepare Standards.

**Caution:** The Human Osteonectin Calibrator is a human source material (isolated from thrombin-activated human platelets) and should be handled as a Biosafety Level 2 material.

MSD supplies a multi-analyte calibrator blend that contains 0.2, 2.0, and 0.2 µg/mL of OCL, ONN, and OPN, respectively. These concentrations reflect the top of the standard curve and this calibrator blend provided should be used as the highest calibrator; no dilution is necessary.

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

- 1) Vortex the calibrator blend provided, briefly, and use it as the highest standard (STD-1).
- 2) Prepare the next standard (STD-2) by transferring 15 µL of the highest standard to 135 µL of Diluent 7. Mix well by vortexing. Repeat 10-fold serial dilutions 5 additional times to generate 7 standards.
- 3) Use Diluent 7 as the zero standard.

### Notes:

*Blocker A solution may be stored at 2–8°C for up to 5 weeks.*

*Detection antibody solution is stable on ice for a few hours.*

*The calibrator dilutions should be prepared immediately before use and kept on ice until use.*



This yields the following calibrator concentrations:

Calibrator	OCL (pg/mL)	ONN (pg/mL)	OPN (pg/mL)
STD-1	200,000	2,000,000	200,000
STD-2	20,000	200,000	20,000
STD-3	2,000	20,000	2,000
STD-4	200	2,000	200
STD-5	20	200	20
STD-6	2.0	20	2.0
STD-7	0.2	2.0	0.2
STD-8	0	0	0

**Notes:**

*Remaining undiluted combined Calibrator stock solutions may be refrozen on dry ice in single use aliquots and stored at  $\leq -70^{\circ}\text{C}$ .*

**Prepare Samples.** Dilute samples with Diluent 7. For serum and plasma samples, MSD recommends a 20-fold sample dilution; however, depending on the sample set under investigation, you may need to use a higher dilution factor. For example, to dilute 20-fold, add 10  $\mu\text{L}$  of sample to 190  $\mu\text{L}$  of Diluent 7.

### Assay Protocol

Begin with a MULTI-SPOT 96-well 7-Spot Human Bone Panel II Plate. No pre-treatment is necessary.

- 1) Add 150  $\mu\text{L}$  of Blocker A solution to each well of the Human Bone Panel II plate. Seal the plate, and incubate for 30 minutes to 1 hour at room temperature.  
Prepare samples and calibrators during this time.
- 2) Wash plates 3 times with 200  $\mu\text{L}$  per well 1X PBS. Add 25  $\mu\text{L}$  of Diluent 7 to all wells. Next, add 25  $\mu\text{L}$  of calibrator dilutions to appropriate calibrator wells, and 25  $\mu\text{L}$  of diluted serum/plasma samples to sample wells. Seal the plate and incubate with shaking for two hours at room temperature.
- 3) Wash plates 3 times with 200  $\mu\text{L}$  per well 1X PBS. Add 25  $\mu\text{L}$  of detection antibody solution to each well. Seal the plate and incubate with shaking for 1 hour at room temperature.
- 4) Wash plates 3 times with 200  $\mu\text{L}$  per well 1X PBS. Add 150  $\mu\text{L}$  of 1X MSD Read buffer T solution to each well and read immediately on the MSD instrument.

*Shaking the plate accelerates analyte capture.*

*Bubbles in the read buffer will interfere with reliable imaging of the plate if carried into the wells.*

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, MSD GOLD, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, S-PLEX, U-PLEX, 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), S-PLEX (design), U-PLEX (design), V-PLEX (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC. © 2015 Meso Scale Diagnostics, LLC. All rights reserved.

