

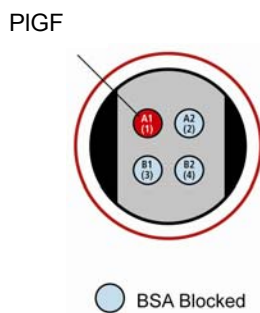
# MSD<sup>®</sup> 96-Well MULTI-ARRAY<sup>®</sup> Human PIGF Assay

The following assay protocol has been optimized for analysis of placental growth factor (PIGF) in human serum and plasma samples.

Storage

## MSD Materials

<input type="checkbox"/> Read Buffer T (4X)	RT
<input type="checkbox"/> Blocker C	2-8 °C
<input type="checkbox"/> MULTI-SPOT <sup>®</sup> 96-well 4 Spot Human PIGF Plate	2-8 °C
<input type="checkbox"/> SULFO-TAG <sup>™</sup> Anti-hPIGF Antibody (100X) <sup>1</sup>	2-8 °C
<input type="checkbox"/> Diluent 7	≤-10 °C
<input type="checkbox"/> Diluent 8	≤-10 °C
<input type="checkbox"/> Diluent 9	≤-10 °C
<input type="checkbox"/> Human PIGF Calibrator (1 µg/mL)	≤-70 °C



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> Some SULFO-TAG labeled detection antibodies may be light-sensitive, so they should be stored in the dark.

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



## Notes:

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

- ❑ Deionized water for diluting Read Buffer
- ❑ Phosphate buffered saline + 0.05% Tween-20 (PBS-T) for plate washing
- ❑ Adhesive plate seals
- ❑ Microtiter plate shaker
- ❑ Plate washer or other efficient multi-channel pipetting equipment for washing 96-well plates
- ❑ Appropriate liquid handling equipment for desired throughput that must accurately dispense 25 and 150  $\mu\text{L}$  into a 96-well microplate

*Read the entire detailed instructions before beginning work.*

### *Protocol at a Glance*

**The protocol can be completed in approximately 5.5 hours** if each reagent is prepared during the preceding incubation. This time can be reduced to **4.5 hours if the blocking reagent is added the night before.**

*The assay protocol was optimized for serum and plasma samples. For significantly different sample matrices, it is recommended to use a calibrator diluent that is similar to the sample matrix (e.g. lysis buffer + carrier protein).*

- Step 1.** Add Blocking Solution, incubate 1-2 hour, wash. (alternatively, block plates overnight at 4 °C).
- Step 2.** Add 25  $\mu\text{L}$  of Diluent 7.  
Add 25  $\mu\text{L}$  of Samples or Calibrator, incubate 2 hours, wash.
- Step 3.** Add 25  $\mu\text{L}$  of Detection Antibody, incubate 2 hours, wash.
- Step 4.** Add 150  $\mu\text{L}$  of Read Buffer, read plate and analyze data.

### *Preparation Instructions*

#### **Thaw Diluents**

Thaw Diluent 7, Diluent 8, and Diluent 9. Vortex briefly. If there is a precipitate, mix gently and warm to room temperature to dissolve. Diluents are stable at 4 °C for one week.

#### **Prepare Calibrator dilutions:**

1. Determine how many Calibrator levels and replicates will be tested in the experiment. Each well will require 25  $\mu\text{L}$  of Calibrator. Thaw Diluent 9 and one vial of Human PIGF Calibrator (1  $\mu\text{g}/\text{mL}$ ). Vortex briefly. Prepare the required Calibrator dilution series using Diluent 9.



**Notes:**

2. A recommended Calibrator dilution procedure is listed below for 4 replicates of 7 Calibrator concentrations spanning a wide range, plus 1 zero-Calibrator point.
  - i. Prepare 1000  $\mu\text{L}$  of PIGF Calibrator at a concentration of 10000  $\text{pg}/\text{mL}$  by adding 10  $\mu\text{L}$  of the PIGF Calibrator at 1  $\mu\text{g}/\text{mL}$  to 990  $\mu\text{L}$  Diluent 9.
  - ii. Prepare 6 additional 1:4 serial dilutions, beginning with the PIGF Calibrator at 10 000  $\text{pg}/\text{mL}$ , by adding 50  $\mu\text{L}$  of the Calibrator to 150  $\mu\text{L}$  Diluent 9.
  - iii. This will create 7 Calibrators with 10000, 2500, 625, 156, 39, 9.8 and 2.4  $\text{pg}/\text{mL}$ . The recommended 8<sup>th</sup> dilution is Diluent 9 alone (e.g. zero Calibrator).
3. Calibrators are stable at room temperature for a few hours. The Calibrator stock solution is stable for one day at 4 °C and for one additional freeze-thaw cycle. Diluent 9 is stable for one week at 4 °C.

**Prepare the 1X Detection Antibody Solution**

- a) In a 15 mL tube combine:
  - 30  $\mu\text{L}$  of 100X SULFO-TAG Anti-hPIGF-1 Antibody
  - 2.97 mL of Diluent 8
- b) This will yield 3 mL of diluted Detection Antibody Solution at the working concentration with sufficient volume for one plate.

*Detection Antibody Solution is stable at RT for a few hours.*

**Dilute Read Buffer:**

- In a 50 mL tube combine (per plate):
- 5 mL 4X Read Buffer T
  - 15 mL deionized water

*Diluted Read Buffer may be kept in a tightly sealed container at room temperature for later use.*



**Notes:**

## *Assay Protocol*

Begin with a MULTI-SPOT 96-well 4 Spot Human PIGF plate.  
No pre-treatment is necessary.

1. Add 150  $\mu$ L/well of Blocker C and incubate at room temperature for 1-2 hours or overnight at 4 °C.
2. Wash plates 3 times with phosphate buffered saline + 0.05%Tween-20 (PBS-T).
3. Add 25  $\mu$ L/well of Diluent 7.
4. Add 25  $\mu$ L/well Calibrator or sample and incubate at room temperature with shaking for 2 hours.
5. Wash plates 3 times with PBS-T.
6. Add 25  $\mu$ L/well of 1X Detection Antibody Solution and incubate at room temperature with shaking for 2 hours.
7. Wash plates 3 times with PBS-T.
8. Prepare SECTOR Imager so that plate can be read immediately after Read Buffer addition.
9. Add 150  $\mu$ L/well 1X Read Buffer T.
10. Analyze immediately with SECTOR Imager.

*Plates may also be blocked overnight at 4°C and stored for up to a week with Blocker C.*

*Avoid bubbles while adding the Read Buffer; it will interfere with accurate reading of the plate.*