

MSD[®] Human Fertility Panel

Application

For the quantitative measurement of LH, FSH, and Progesterone in human serum

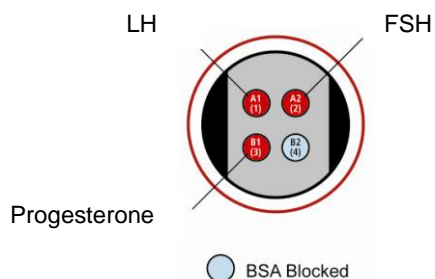
Storage

MSD Materials

<input type="checkbox"/> Read Buffer T (4X)	RT
<input type="checkbox"/> MULTI-SPOT [®] 96-well Human Fertility Panel Plate	2-8 °C
<input type="checkbox"/> SULFO-TAG [™] LH Detection Antibody (500 µg/mL)	2-8 °C
<input type="checkbox"/> SULFO-TAG FSH Detection Antibody (500 µg/mL)	2-8 °C
<input type="checkbox"/> SULFO-TAG Progesterone (1 µg/mL)	2-8 °C
<input type="checkbox"/> Diluent 22	≤-10 °C
<input type="checkbox"/> Fertility Calibrator Set ¹ (combined set of 8 concentrations, 125 µL each)	-20 °C

Other Materials & Equipment (not supplied)

- Deionized water for diluting Read Buffer
- Phosphate Buffered Saline for plate washing
- Automated 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 µL into a 96-well micro plate



The SECTOR[®] Imager data file will identify spots according to their well location, not by the coated capture antibody name.

¹ The Calibrator blend in this kit is derived from human source material which has been tested and found to be negative for HIV-1 and 2, Hepatitis B, and Hepatitis C. 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 OR THERAPEUTIC PROCEDURES.



Protocol at a Glance

Notes:

The protocol can be completed in approximately 1.5 hours.

Read the entire detailed instructions before beginning work.

1. Add Detection Antibody Reagent to each well.
2. Add Calibrator or sample and incubate for 1 hour with shaking.
3. Wash.
4. Add Read Buffer and Read plate.

Preparation Instructions

Prepare Calibrators:

1. Thaw one set of Calibrators: each set contains a sufficient volume for four replicates of each Calibrator (25 μ L/well, 125 μ L provided). Once thawed, Calibrators should be refrigerated (for up to 8 hours) if not used immediately. Calibrators can tolerate one additional freeze/thaw cycle.
2. The Calibrators contain all three fertility markers: the concentration of each marker is listed on the Calibrator tube label (units: mIU/mL for LH and FSH; ng/mL for progesterone). At least two replicates of each Calibrator should be included on each plate to produce a well-defined calibration curve.

Prepare Detection Antibody Reagent:

1. Determine total number of wells in the experiment. Each well will require 25 μ L of Diluent 22. Thaw an appropriate amount of Diluent 22 for the experiment. Diluent 22 can be refrozen up to 3 times.
2. Prepare Detection Antibody Reagent by diluting all three SULFO-TAG reagents **together into a single volume** of Diluent 22 as follows:
 - Dilute LH Antibody (stock – 500 μ g/mL) to a final concentration of 2 μ g/mL
 - Dilute FSH Antibody (stock – 500 μ g/mL) to a final concentration of 12 μ g/mL
 - Dilute labeled Progesterone (stock – 1 μ g/mL) to a final concentration of 2 ng/mL
3. Detection Antibody Reagent is stable at room temperature for a few hours.



Notes:

Dilute Read Buffer:

1. Determine total number of wells in the experiment. Each well will receive 150 μ L of Read Buffer T. Prepare an extra 20%.
2. Dilute 4X Read Buffer T to 1X with deionized water.
3. Diluted Read Buffer may be stored at room temperature for later use.

Assay Protocol

Begin with a MULTI-SPOT 96-well Human Fertility Panel Plate.

1. Add 25 μ L/well of Detection Antibody Reagent (solution of Diluent 22 containing two diluted SULFO-TAG antibodies and labeled progesterone).
2. Add 25 μ L/well Calibrator or sample and incubate at room temperature with shaking for 1 hour.
3. Prepare SECTOR instrument such that the plate can be read immediately following Read Buffer addition.
4. Wash plates 3 times with PBS.
5. Add 150 μ L/well 1X Read Buffer T. Avoid bubbles. The use of an electronic multi-pipettor at moderate speed setting is recommended.

Note that bubbles introduced to the well during Read Buffer addition will interfere with reliable imaging of the plate.

