

R-PLEX™ Antibody Set

Human FAP- α /SEPR

Cat. No: F21YI-3, 5-plate size

F21YI-8, 50-plate size

Ordering Information

MSD Customer Service
Phone: 1-240-314-2795
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com
www.mesoscale.com/support

Scientific Support

Phone: 1-240-314-2798
Email: ScientificSupport@mesoscale.com
www.mesoscale.com/support

Company Address

MESO SCALE DISCOVERY®
A division of
Meso Scale Diagnostics, LLC.
1601 Research Boulevard
Rockville, MD 20850-3173 USA

www.mesoscale.com®

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

Contents and Storage

Description

| Description | No. of vials | Vol. /vial | Storage |
|---|--------------|---------------|----------------------------|
| Human FAP- α /SEPR Calibrator, 20X | 5 | 50 μ L | $\leq -70^{\circ}\text{C}$ |
| Biotin Human FAP- α /SEPR Capture Antibody | 1 | 1,250 μ L | 2-8 $^{\circ}\text{C}$ |
| SULFO-TAG™ Human FAP- α /SEPR Detection Antibody, 100X | 1 | 375 μ L | 2-8 $^{\circ}\text{C}$ |

Note: Calibrators and antibodies are shipped separately.

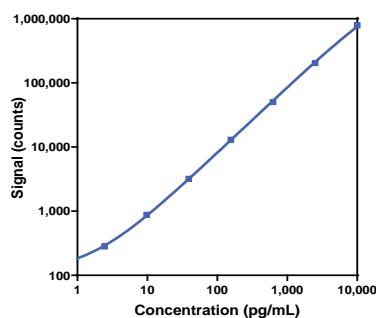
R-PLEX Antibody Sets provide an expanded menu of electrochemiluminescence assays for biomarker discovery and development. The sets include a matched biotinylated capture and SULFO-TAG conjugated detection antibody pair and calibrator for the quick and easy development of highly sensitive immunoassays on a variety of MSD® plates including U-PLEX® MULTI-SPOT plates and MSD GOLD™ Small Spot Streptavidin plates.

Fibroblast Activation Protein- α (FAP- α , seprase) is a member of the serine protease family and is a type 2 membrane-bound glycoprotein. It has similar activity as DPPIV peptidase and collagenases. FAP- α has been associated with enhancing tumor activity in cell lines. Elevated levels of FAP- α in the stroma indicates poor prognosis in epithelial cancer. Sites of tissue remodeling generally have high FAP- α .

Assay Performance

R-PLEX Antibody Sets can be easily adapted to develop singleplex and multiplex assays on the MSD platform. MSD GOLD Small Spot Streptavidin plates can be used for singleplex assay development, whereas multiplex assay formats can be performed using U-PLEX MULTI-SPOT® plates.

Representative data from the R-PLEX Human FAP- α /SEPR Antibody Set tested on MSD's U-PLEX platform using Diluent 101 as assay diluent and Diluent 37 as antibody diluent is presented below. The lower limit of detection (LLOD) is a calculated concentration corresponding to the signal 2.5 standard deviations above the background (zero calibrator). The performance of your assay may vary depending on the diluent used, the sample type, and the other analytes when testing in a multiplexed format.



| Top-of-Curve Conc. (pg/mL) | LLOD (pg/mL) |
|----------------------------|--------------|
| 10,000 | 1.0 |

For developing individual assays using R-PLEX Antibody Sets, please refer to the R-PLEX Antibody Sets Singleplex Assays product insert. For multiplex assay development using R-PLEX Antibody Sets or a combination of R-PLEX and U-PLEX Antibody Sets, refer to the U-PLEX Development Pack product insert. The detailed instruction manuals as well as the R-PLEX assay datasheets are available at www.mesoscale.com/R-PLEX-documents.

Sample Dilution

The R-PLEX Human FAP- α /SEPR assay has been tested with the following sample type(s) at the specified dilution in assay diluent.

| Sample Type | Recommended Fold Dilution |
|--------------|---------------------------|
| Human Plasma | 100 |

The assay may be used for other sample types. We recommend performing a dilution series and testing for diluent compatibility when measuring a different sample type.

R-PLEX Antibody Set

Source

| | |
|------------------|--|
| Calibrator | Recombinant Fibroblast Activation Protein- α |
| Antibodies | Capture antibody - mouse monoclonal, detection antibody - sheep polyclonal |
| Assay generation | A |

Recommended Assay Components

The following items are recommended for developing assays using R-PLEX Antibody Sets.

Plates for Singleplex Assay Development: All plates are available in 1, 5, 30, 120, and 510 plate sizes.

| Description | Catalog No. |
|---|-------------|
| MSD GOLD 96-well Small Spot Streptavidin Plates | L45SA |

Plates for Multiplex Assay Development: The U-PLEX platform allows you to create your own multiplex assays. Using two simple tools, a 10-spot U-PLEX plate and unique Linkers, you can build custom multiplex panels for any combination of analytes. U-PLEX Development Pack includes U-PLEX Plates, Linkers, Stop Solution, and Read Buffer T. All development packs are available in 1, 5, and 25-plate pack sizes. You can choose the appropriate Development Pack according to the number of assays you plan to multiplex, e.g., use a U-PLEX 4-Assay Development Pack if you are developing a 4-plex assay.

| Description | Catalog No. |
|---------------------------------|-------------|
| U-PLEX 2-Assay Development Pack | K15227N |
| U-PLEX 3-Assay Development Pack | K15228N |
| U-PLEX 4-Assay Development Pack | K15229N |
| U-PLEX 5-Assay Development Pack | K15230N |
| U-PLEX 6-Assay Development Pack | K15231N |

| Description | Catalog No. |
|----------------------------------|-------------|
| U-PLEX 7-Assay Development Pack | K15232N |
| U-PLEX 8-Assay Development Pack | K15233N |
| U-PLEX 9-Assay Development Pack | K15234N |
| U-PLEX 10-Assay Development Pack | K15235N |

Note: R-PLEX assays can be combined with U-PLEX assays to build multiplex panels. View the R-PLEX Assay Designer at www.mesoscale.com/R-PLEX to customize your multiplex panel.

Buffers:

| Description | Catalog No. |
|-------------------------------|-------------|
| MSD GOLD Read Buffer (200 mL) | R92TG-1 |
| MSD Wash Buffer (20X, 100 mL) | R61AA-1 |

Diluent Compatibility

The R-PLEX Human FAP- α /SEPR Antibody Set has been tested with the following diluents.

| Diluent Type | Diluent Name | Catalog No. | Size |
|------------------|--------------|-------------|--------|
| Assay Diluent | Diluent 101 | R51AD-3 | 50 mL |
| | | R51AD-7 | 700 mL |
| Antibody Diluent | Diluent 37 | R50AF-3 | 25 mL |
| | | R50AF-6 | 125 mL |

To run five plates, 50 mL of assay diluent and 40 mL of antibody diluent are required. Additional assay diluent may be required for sample dilution. Testing of different diluents can help optimize assays for specific experimental conditions. MSD offers a range of assay and antibody diluents for separate purchase. Please view our website at www.mesoscale.com/R-PLEX for more details on compatible diluents for R-PLEX assays.

The R-PLEX Human FAP- α /SEPR Antibody Set is for research use only.