

# Immobilization of Non-Adherent Cells with Cell-Tak for Assay on the Agilent Seahorse XFe/XF96 or XFp Analyzer

## Technical Overview

### Introduction

Cell-Tak cell and tissue adhesive may be used to prepare adherent monolayer cultures of biological samples normally grown in suspension such as lymphocytes and platelets, for assay on the Agilent Seahorse XFe/XF96 or Agilent Seahorse XFp Analyzer<sup>1-7</sup>. Cell-Tak is a nonimmunogenic extracellular matrix protein preparation isolated from the marine mussel, *Mytilus edulis*<sup>8</sup>.



The following protocol (Figure 1) is for cells grown in suspension that do not naturally settle to the bottom of the microplate well under gravity, thus requiring centrifugation to settle down. For cells that naturally settle down to the bottom of the well, skip to Seeding Cells in Cell-Tak-Coated Plates step 11.

## Materials

- Agilent Seahorse XF Base Medium (p/n 102353-100 for 2 L; 103193-100 for 100 mL)
- Agilent Seahorse XF96 Cell Culture Microplate (p/n 101085-004), or Agilent Seahorse XFp Cell Culture Miniplate (p/n 103025-100)
- Corning Cell-Tak Cell and Tissue Adhesive, 1 mg (Corning, Cat. # 354240)
- Sodium bicarbonate (NaHCO<sub>3</sub>) (Sigma, Cat. # S5761)
- Sodium hydroxide (Sigma, Cat # 38215)
- Tissue Culture Grade Sterile Water (Invitrogen, Cat. # 15230)
- Water bath set at 37 °C
- Pipettors (single or multichannel)

- Optional, depending on cell type:** Benchtop centrifuge with swing-bucket rotor equipped with plate carriers. Example: Eppendorf Centrifuge 5810R.

## Preparation of Cell-Tak-Coated Plates

Follow the manufacturer's Basic Absorption Coating Protocol, and refer to the Coating Procedure for Multiple Well Plates outlined in the Instructions for Use<sup>8</sup>. Agilent has identified the following reference points and exceptions helpful to adapt the Coating Procedure for use with Agilent Seahorse XF96 Cell Culture Microplates and Seahorse XFp Cell Culture Miniplates:

- The optimal Cell-Tak solution concentration for Agilent Seahorse XF96 Cell Culture Microplate or Agilent Seahorse XFp Cell Culture Miniplate is 22.4 µg/mL.
- Prepare 2.5 mL of this solution for the Seahorse XFe/XF96, or 0.25 mL of this solution for the Seahorse XFp Analyzer. Refer to the Manufacturer's protocol to prepare this solution.
- Apply 25 µL of the solution to each well for 20 minutes at room temperature.

- Wash each well twice using 200 µL of sterile water.
- Cell-Tak-coated Seahorse XF96 Cell Culture Microplates and Seahorse XFp Cell Culture Miniplates may be stored for up to 1 week at 4 °C.
- Cell-Tak coated XF96 Cell Culture Microplates and XFp Cell Culture Miniplates must be allowed to warm to room temperature in the hood before cell seeding.

**Note:** Per manufacturer's Instructions for Use, do not pre-incubate serum-containing medium in the Cell-Tak-coated wells prior to cell seeding, as this may result in a loss of adhesion.

## Seeding Cells in Cell-Tak-Coated Plates

**Note:** Optimal cell density may vary between cell types. Agilent recommends optimizing cell density parameters prior to beginning the assay to ensure reproducible results.

The following protocol (Figure 1) describes seeding one Seahorse XF96 Cell Culture Microplate or Seahorse XFp Cell Culture Miniplate. To balance the centrifuge, create a dummy plate by adding 50 µL of water to each well.

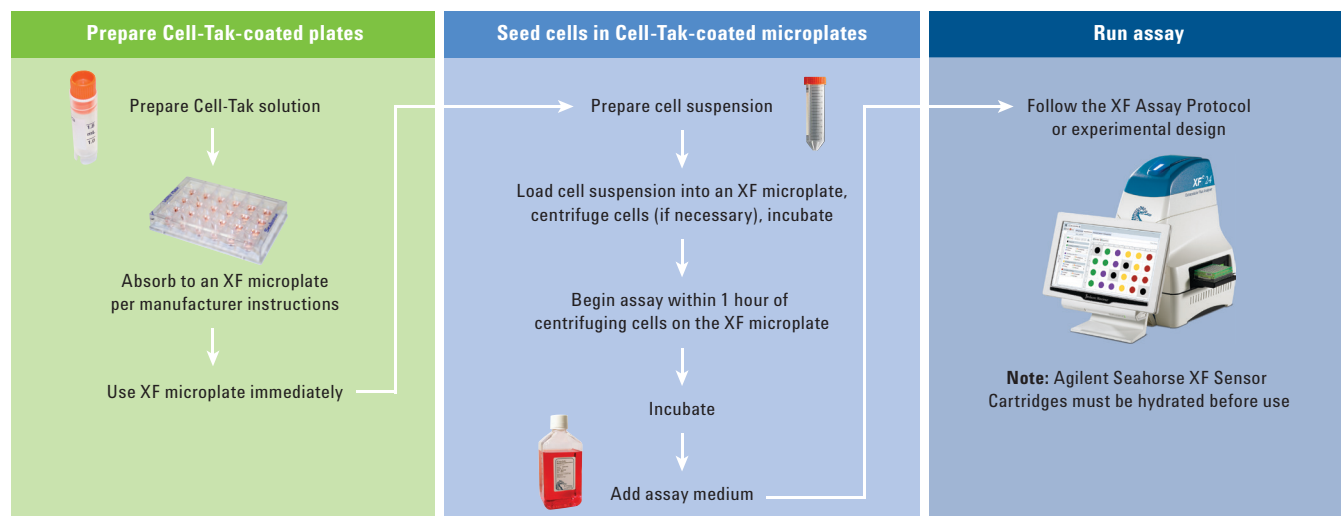


Figure 1. Protocol flow chart.

1. Prepare assay medium, by supplementing Seahorse XF Base Medium as required by the experimental conditions. Warm medium in a 37 °C water bath.
2. For one Seahorse XF96 Cell Culture Microplate or Seahorse XFp Cell Culture Miniplate, transfer an appropriate volume of cell suspension from the growth vessel to a conical tube. To calculate the total number of cells needed, multiply the desired number of cells per well times 100 wells or 10 wells for the Seahorse XFe/XF96 or XFp, respectively. (For example, 150,000 cells per well  $\times$  100 wells =  $1.5 \times 10^7$  cells).
3. Centrifuge cells at room temperature at  $200 \times g$  for 5 minutes.
4. While cells are being centrifuged, pipette 50  $\mu$ L assay medium into background/control wells of the room-temperature Cell-Tak-coated Seahorse XF96 Cell Culture Microplates or Seahorse XFp Cell Culture Miniplates.
5. Remove supernatant from the centrifuged conical tube.
6. Resuspend cells in warmed assay medium to the desired concentration of cells per well in 50  $\mu$ L of assay medium. (For example,  $1.5 \times 10^5$  cells per well is desired, resuspend cells in a volume that results in  $1.5 \times 10^5$  cell/50  $\mu$ L or  $3.0 \times 10^6$  cells/mL).
7. Change centrifuge settings to zero braking.
8. Transfer the cell suspension to a sterile tissue culture reservoir, or pipette from the conical tube.
9. Pipette 50  $\mu$ L of the cell suspension along the side of each well, except for background/control wells. Agilent recommends using a multichannel pipette.
10. Centrifuge the cells at  $200 \times g$  (zero braking) for 1 minute. Ensure that the centrifuge is properly balanced. For XFp Analyzer users, Agilent recommends using the Agilent Seahorse XFp Carrier Tray to centrifuge the Seahorse XFp Cell Culture Miniplates. For more details, refer to the Basic Procedure: Seeding Suspension Cells in XFp Cell Culture Miniplates.
11. Transfer plates to a 37 °C incubator not supplemented with CO<sub>2</sub> for 25–30 minutes to ensure that the cells have completely attached. Visually confirm that most of the cells are stably adhered to the culture surface.
12. Slowly and gently, add 130  $\mu$ L warm assay medium along the side of each well. Take care to avoid disturbing the cells.
13. Observe the cells under the microscope to check that cells are not detached.
14. Return the cell plates to the incubator for 15–25 minutes.
15. After 15–25 minutes, the cell plates are ready for assay. Total time following centrifugation should be no greater than 1 hour for best results.
16. Place the cell plate in either the Seahorse XFe/XF or XFp Analyzer, following calibration.
17. Proceed following the assay protocol.

## Notes

This protocol specifies the full-plate seeding of a single Seahorse XF96 Cell Culture Microplate or Seahorse XFp Cell Culture Miniplate. If more than one plate is desired, increase the volumes and total cell numbers required proportionately. Agilent recommends seeding two plates when beginning work with a cell line for additional practice with step 12 (addition of medium without disrupting cells).

This cell seeding protocol was developed by Agilent Seahorse Bioscience scientists, and is applicable exclusively to cells cultured in the Seahorse XF96 Cell Culture Microplates or Seahorse XFp Cell Culture Miniplates coated with Cell-Tak, and is intended for analysis using a Seahorse XFe/XF96 or XFp Analyzer.

## References

1. Wang, R.; *et al. Immunity*. **2011**, *35*(6), 871-82 (T-lymphocytes).
2. Capasso, M.; *et al. Nat. Immunol.* **2010**, *11*(3), 265-72 (B-lymphocytes).
3. Avila, C.; *et al. Exp. Clin. Endocrinol. Diabetes*. **2011**, DOI: 10.1055/s-0031-1285833 2011 (platelets).
4. Stackley, K.; *et al. PLoS One* **2011**, *6*(9), e25652 (zebrafish embryos).
5. Rogers, G.; *et al. PLoS One* **2011**, *6*(7), e21746 (isolated mitochondria).
6. Bulua, A. C.; *et al. J. Exp. Med.* **2011**, *208*(3), 519-33 (PBMC).
7. Saha, A.; *et al. Gut* **2010**, *59*(7), 874-881 (AGS adenocarcinoma cells)
8. Corning Cell-Tak Cell and Tissue Adhesive Instructions for Use.

## Further Reading

Agilent Seahorse Bioscience Application Note (available online):

*Bioenergetic analysis of suspension cells: hematopoietic stem cells and lymphocytes.* A real-time assay that quantifies the ATP and biosynthetic demands of immune cell proliferation, differentiation, and effector function.

[www.agilent.com](http://www.agilent.com)

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

This information is subject to change without notice.

© Agilent Technologies, Inc., 2016  
Published in the USA, December 1, 2016  
5991-7153EN



**Agilent Technologies**