

Washing Adherent Cells in Agilent Seahorse XF24 Cell Culture Microplates

Basic Procedure

Before performing an Agilent Seahorse XF Assay, growth medium must be replaced with a suitable assay medium (generally, medium without bicarbonate buffer or serum and low/no phenol red content). This procedure describes replacing the growth medium with assay medium for adherent cells grown in XF24 Cell Culture Microplates prior to being assayed using an Agilent Seahorse XFe24/XF24 Analyzer.

- Warm the assay medium to 37°C.
 For choosing and preparing the appropriate assay medium, please see http://www.agilent.com/cs/library/selectionguide/public/5991-7878EN. pdf and http://www.agilent.com/cs/library/usermanuals/public/XFe24_DAY_OF_MEDIA_PREP.pdf or http://www.agilent.com/cs/library/usermanuals/public/XF24_DAY_OF_MEDIA_PREP.pdf.
- 2. Retrieve the cell culture microplate from the CO_2 incubator.
- 3. Look at cells under the microscope to:
 - a. Confirm cell health, morphology, seeding uniformity and purity (no contamination).
 - b. Ensure cells are adhered, and no gaps are present.
 - c. Make sure no cells were plated in the background correction wells
- 4. Wash cells with assay medium
 - a. Remove all but 50 µL of the culture medium from each well.
 - b. Rinse cells two times with 1 mL of assay medium, leaving behind $50 \mu l$ after each wash.
 - c. Add 450 μ L of assay medium to each well for a final volume of 500 μ L/well.
- 5. Look at cells under the microscope to ensure that cells were not disturbed or washed away.
- 6. Place the plate in a 37°C incubator without CO₂ for 45-60 minutes prior to the assay.

NOTE: Incubating the cell plates without CO₂ allows outgassing from the plate and is required for accurate ECAR measurements.



www.agilent.com/en-us/promotions/seahorse-xf-technology

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

Product specifications and descriptions in this document are subject to change without notice.

© Agilent Technologies, Inc., 2014, 2017 Published in the USA, June 30, 2017

