



# Agilent Seahorse XF<sup>e</sup>96 Analyzer Materials List

Materials required and recommended for a successful XF<sup>e</sup>96 assay.

## Required

The following are the materials required to perform an XF<sup>e</sup>96 assay.

1. Agilent Seahorse XF<sup>e</sup>96 Analyzer with Controller
2. Non-CO<sub>2</sub> Incubator (37°C)
3. Agilent Seahorse XF<sup>e</sup>96 FluxPak (Part # 102416-001) or Agilent Seahorse XF<sup>e</sup>96 FluxPak mini (Part # 102601-100)
4. Calibrated pH Meter
5. Inverted Phase Contrast Microscope
6. P20 and P200 Pipettes and tips
7. Appropriate XF Assay Media (for guidance in choosing and preparing the proper assay media, see: <http://www.agilent.com/cs/library/selectionguide/public/5991-7878EN.pdf> and [http://www.agilent.com/cs/library/usermanuals/public/XFe96\\_DAY\\_OF\\_MEDIA\\_PREP.pdf](http://www.agilent.com/cs/library/usermanuals/public/XFe96_DAY_OF_MEDIA_PREP.pdf), respectively).



## Recommended

The following are materials recommended to help achieve optimal results.

1. Uninterrupted power source
2. Internet access
3. Balance
4. Stir plate
5. Cell counter/Hemocytometer
6. Water bath set at 37°C
7. Microcentrifuge
8. Touch vortex
9. Centrifuge with adaptors for cell culture plates
10. Reservoirs
11. 15 and 50 mL conical tubes
12. Sterile filter bottles (0.22 µm filter) and cap

## Available Kits and Reagents

A variety of materials is offered by Agilent Technologies to assist with designing assays to examine cellular metabolism, including assay kits, reagents, media and special assay plate designs.

If you have any questions about which kits, reagents and plates may be appropriate for your XF experiments, please visit the Agilent website: [http://www.agilent.com/en-us/products/cell-analysis-\(seahorse\)/seahorse-xfe-consumables](http://www.agilent.com/en-us/products/cell-analysis-(seahorse)/seahorse-xfe-consumables), or contact technical support: [seahorse.support@agilent.com](mailto:seahorse.support@agilent.com).

Learn more  
[www.agilent.com/en-us/promotions/seahorse-xf-technology](http://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



**Agilent Technologies**