

## Recommendations for Decontamination of an Agilent Seahorse XF Analyzer



The use of vaporized hydrogen peroxide (VHP) for decontamination of Agilent Seahorse XF instruments, including XF, XFe and XFp platforms, has been validated. The use of VHP will not result in any damage to the instrument nor void the instrument warranty if performed as recommended below.

It is recommended that the decontamination process be carried out by a certified vendor. For specific vendor recommendations, please email Cell Analysis technical support at [cellanalysis.support@agilent.com](mailto:cellanalysis.support@agilent.com).

Decontamination is carried out according to the vendor's standard protocol. In brief, the analyzer is completely covered and taped with a plastic bag with the instrument's cover(s) off/opened. VHP is injected into the sealed bag at controlled humidity ( $RH \leq 60\%$ ) and air flow for 1 hour. The bag is then opened and the vapor allowed to dissipate.

Pre- and post-decontamination XF assays have been performed by Agilent to insure proper instrument performance. VHP exposure does not cause any physical damage to the Seahorse XF Analyzer components nor the functional aspects of the instrument when performed as recommended above.



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

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

This information is subject to change without notice.