

Xuri™ W25

Addendum

About this addendum

This addendum describes changes to existing information. This document is an addendum to the documents listed below.

- *Xuri™ Cell Expansion System W25 Operating Instructions 29064612*
- *Xuri Cell Expansion System W25 User Manual 29064622*

Change of product name

The product name Xuri Cell Expansion System W25 has been changed to Xuri W25.

The change of product name is not reflected in the Operating Instructions and User Manual. Hence, the product name on the system labels, instrument labels, and packaging labels does not align with the product name in the *Xuri Cell Expansion System W25 Operating Instructions 29064612* and *Xuri Cell Expansion System W25 User Manual 29064622*.

The Operating Instructions and User Manual are valid for Xuri W25.



cytiva.com

Cytiva and the Drop logo are trademarks of Life Sciences IP Holdings Corp. or an affiliate doing business as Cytiva.

Xuri is a trademark of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

All other third-party trademarks are the property of their respective owners.

© 2022 Cytiva

For local office contact information, visit cytiva.com/contact

29714514 AA V:1 01/2022