

Sartorius Stedim Biotech GmbH 37070 Goettingen Germany

For additional information, please contact your local Sartorius Stedim Biotech representative.i.V.

Sartorius Stedim Biotech GmbH
August-Spindler-Straße 11
37079 Göttingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius-stedim.com

April 1st 2013

Change Notification UniVessel® SU

Introduction

As an international leading supplier of process technologies to the bio-pharmaceutical industry, Sartorius Stedim Biotech is committed to a continuous improvement of its products and services. A defined change control procedure is followed whenever changes to existing products are implemented.

This change notification serves to notify all users of the UniVessel® SU product family of changes to the optical UniVessel® SU vessel holder and the UniVessel® SU 2L culture vessels.

Description of the Changes

In order to harmonize our single-use sensor technology platform for bioreactors from process development to production, we have developed a new opto-electronic device for non-invasive measurement of pH and DO in the UniVessel® SU product family. This new opto-electronic device is using the same measurement technology as already established in the BIOSTAT® RM and BIOSTAT® STR product family.

The product changes include a UniVessel® SU vessel holder with the new integrated opto-electronics and new single-use sensor patches on the UniVessel® SU 2L culture vessels. For both products new article numbers have been implemented.

Affected Products

The current optical UniVessel SU vessel holder (DZ-----VHO) will be replaced by the new optical UniVessel® SU vessel holder identified with order number DZ-----VHC.


The UniVessel® SU 2L (CC1) culture vessels (DU0002LL-SS----V) will be replaced by UniVessel® SU 2L (CC2) culture vessels identified with order number DU0002LL-SS01--V with the new single-use sensor patches.

Please note that the new opto-electronics in the UniVessel® SU vessel holder (DZ-----VHC) will only work in combination with the new UniVessel® SU 2L (CC2) culture vessels (DU0002LL-SS01--V). Existing optical UniVessel® SU vessel holders (DZ-----VHO) will be replaced free of charge, users will be contacted by Sartorius Stedim Biotech to ensure a smooth replacement.

Implementation of the Change

Sartorius Stedim Biotech will discontinue production of the old optical UniVessel SU vessel holder (DZ-----VHO) and the UniVessel® SU 2L (CC1) culture vessels (DU0002LL-SS----V) effective immediately.

The current stock of UniVessel® SU 2L (CC1) culture vessels (DU0002LL-SS----V) can be ordered for use with classical pH and DO sensors. Once consumed, orders will be automatically changed without further notification to the new article UniVessel® SU 2L (CC2) culture vessel (DU0002LL-SS01--V).


Alan Burns
VP Global Quality Systems
Sartorius Corp.


Stefan Schlack
Senior VP Marketing
Sartorius Stedim Biotech GmbH