

Replace Conductivity monitor

Instructions

Scope

Instructions how to replace Conductivity monitor C9 (28956495).

Instruction

The instruction below describes how to replace the conductivity monitor in the instrument.



WARNING

Disconnect power. Always switch off power to the instrument before replacing any component on the instrument, unless stated otherwise in the user documentation.

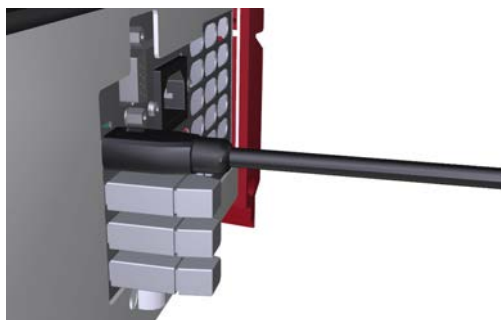
Step	Action
1	Disconnect power from the instrument by switching off the instrument with the Power switch located on the left hand side of the instrument.

Step	Action
------	--------

- | | |
|---|--|
| 2 | Disconnect tubing from the old conductivity monitor. |
|---|--|



- | | |
|---|--|
| 3 | At the back of the instrument, disconnect the conductivity monitor cable . |
|---|--|



- | | |
|---|--|
| 4 | Remove the conductivity monitor from the rail. |
|---|--|

- | | |
|---|--|
| 5 | <ul style="list-style-type: none">• Attach the new monitor to the rail.• Connect the cable to the instrument.• Connect tubing. |
|---|--|

Page intentionally left blank



cytiva.com/akta

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

AKTA™ 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.

© 2020–2021 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

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

28968653 AB V:4 01/2021