

1.17247.0001

# RQcheck

## set for RQflex® 20

For instructions for use, please turn over the page!

Recommended duration for the use of the check strip: 2 years.

Please replace the set immediately if the strip is permanently stained or discolored.

MilliporeSigma is the U.S. and Canada Life Science business of Merck KGaA, Darmstadt, Germany.

© 2025 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. MilliporeSigma, the vibrant M, Supelco, Sigma-Aldrich, and Reflectoquant are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

02/2025



### **Contents:**

**1 grey check strip**

**1 barcode strip "RQcheck 20"**

#### Procedure:

1. Switch on the instrument by pressing the [Power ON/OFF] button.
2. Select [Quality Assurance] - [RQcheck] - [OK].
3. Read the barcode strip into the barcode reader.
4. Now insert the check strip for RQcheck set for RQflex® 20 into the strip adapter (see illustration on front page).
5. Press the [Measure] button.
6. "RQcheck passed!" is shown on the display and the results are automatically stored.
7. Pressing the [OK] button leads forward to [RQcheck Results].

For details see Operating Manual.

#### Trouble shooting:

If an error message is shown on the display and the RQcheck is not passed, clean the strip adapter carefully and recalibrate the instrument using the recalibration set (see RQflex® 20 Operating Manual).

Repeat the check procedure using RQcheck set for RQflex® 20.

If the error message is shown again, please call our technical service department.