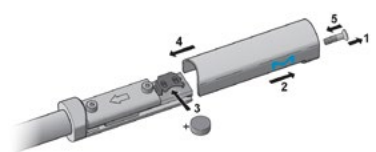


Quick Guide

Luminometer LED Pen Overview

1. Battery replacement



1. Remove the screw
2. Slide the cover
3. Insert the battery ("+" facing up)
4. Close the cover
5. Put back the screw

2. Powering



OFF

ON

3. Battery status



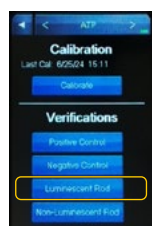
OK

Replace battery*

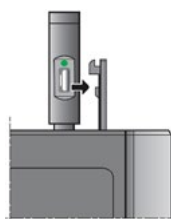
Do not use
Replace battery

*Once the Red Icon shows up, please change the battery.

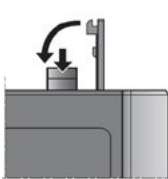
4. Verification mode: Luminescent



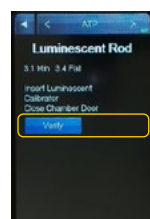
LUMINESCENT



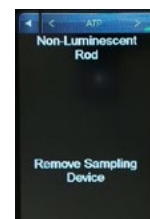
SWITCH ON



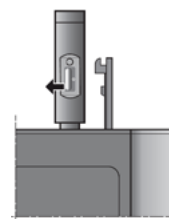
INSERT, CLOSE



SELECT VERIFY

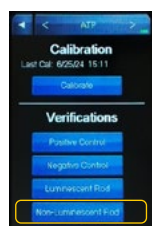


Remove Sampling Device

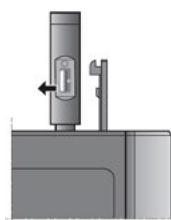


OPEN, REMOVE,
SWITCH OFF

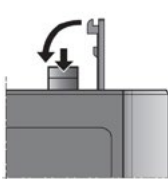
5. Verification mode: Non-Luminescent



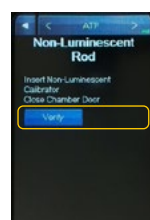
NON-LUMINESCENT



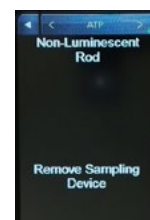
SWITCH OFF



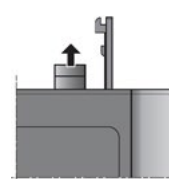
INSERT, CLOSE



SELECT VERIFY



Remove Sampling Device



OPEN, REMOVE

❗ If « **Baseline reading required** » is displayed, remove the LED PEN and follow on-screen prompts.

For full User Guide including usage guidelines, Verification and Calibration steps, cleaning and storage recommendation, visit: www.sigmaaldrich.com



The life science business of Merck KGaA, Darmstadt, Germany operates as MilliporeSigma in the U.S. and Canada.

© 2024 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. The vibrant M, Millipore, Sigma-Aldrich and MVP ICON 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.

VM_UG14066EN | 2024 - 59987 | 12/2024

