

Unveiling the New LED Verification Pen (LTHPEN01)

The newly developed LED Pen enhances the operational efficacy of the MVP ICON system. This innovative product serves as a replacement for the ATP Calibrator Pen (64012) within our Hygiene Monitoring Portfolio.

Advantages of the LED Pen

Calibration and Verification: The LED Pen facilitates both calibration and verification processes for the MVP ICON, ensuring optimal convenience and reliability in performance assessments.

Dual Control Measurement: This single pen is capable of measuring both Negative and Positive Controls, streamlining the verification process.

User-Friendly Design:

Battery Status Indicator: The device features a clear battery status indication system, utilizing green and red signals for easy identification of battery life.

Standard Battery Type: It operates on a standard button battery, allowing for widespread compatibility and ease of replacement.

Battery Replacement Protocol: The design includes a straightforward battery replacement procedure to minimize downtime.

Regulatory Compliance: The LED Pen adheres to international regulatory standards, including CE, EAC, and UK CA certifications, ensuring its suitability for global markets.

Then...



Pen Configuration: One pen is designated for non-luminescent applications, while the other is specifically designed for luminescent light output.

Light Emission Source: Utilizes a low-level C14 source for light emission.

Shelf Life: The product has a shelf life of five years.

Now...



Pen Configuration: A single pen is designed for both non-luminescent and luminescent light output.

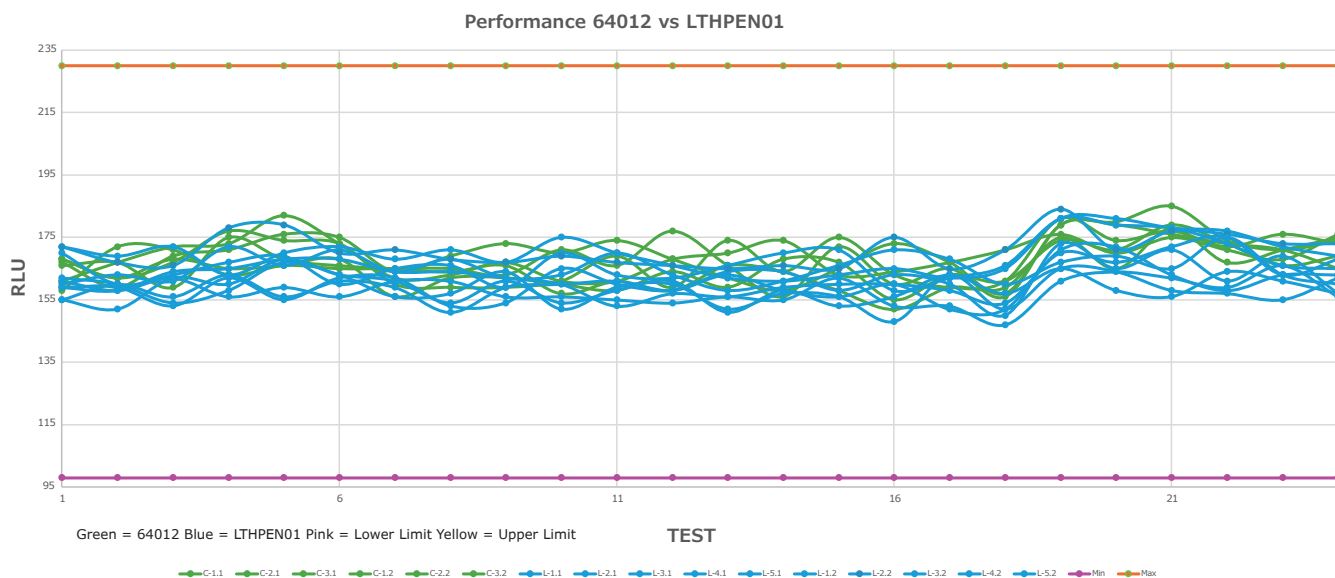
Light Emission Source: Features an electronic-based light emission using LED technology.

Shelf Life: The product has a shelf life of five years.

For more information, please refer to the USER Guide or Quick Guide.



To eliminate variability in the performance of the MVP ICON when calibrated with the new LED Pen, a comprehensive assessment was conducted. This evaluation involved comparing multiple units of the Calibrator Set and the LED Pen across several MVP ICON units.



Calibrator Set (64012), L – LED Pen (LTHPEN01)

The graph clearly demonstrates that the performance of both the Calibrator Set and the new LED Pen falls within the established specified range. Furthermore, the results indicate no significant difference in performance between the two units, as evidenced by the overlapping data obtained from multiple tests conducted.

Need more training? Please register at our LEARN@M platform and get the video training for the LED PEN and other BioMonitoring products.



Ordering Information

LTHPEN01 - LED Verification Pen - 1 unit

To place an order or receive technical assistance

In Europe, please call Customer Service:

France: 0825 045 645 Spain: 901 516 645 Option 1
Germany: 069 86798021 Switzerland: 0848 645 645
Italy: 848 845 645 United Kingdom: 0870 900 4645

For other countries across Europe, please call: +44 (0) 115 943 0840

Or visit: [MerckMillipore.com/offices](https://www.MerckMillipore.com/offices)

For Technical Service visit: [MerckMillipore.com/techservice](https://www.MerckMillipore.com/techservice)

[MerckMillipore.com](https://www.MerckMillipore.com)

We have built a unique collection of life science brands with unrivalled experience in supporting your scientific advancements.

Millipore® Sigma-Aldrich® Supelco® Milli-Q® SAFC® BioReliance®

Merck KGaA
Frankfurter Strasse 250
64293 Darmstadt, Germany

