

Lot specific information

Product: Permafluor E+  
 Lot number: 90-23435

Mfg. Date: 23-OCT-2023  
 Exp. Date: 01-MAY-2026

Information

The product has been analyzed and is certified to have met the specifications.

| Test                              | Temperature | Results                       |
|-----------------------------------|-------------|-------------------------------|
| Appearance                        |             | Clear blue fluorescent liquid |
| Quench parameter                  |             | 78.0%                         |
| Quench parameter with Soluene-350 |             | 56.0%                         |
| Background with Soluene-350       |             | 24 cpm                        |

This Certificate of Analysis has been produced electronically and is valid without a signature.  
 For questions regarding this Certificate of Analysis, please contact your customer service representative.