

Lot specific information

Product: Permafluor E+ Mfg. Date: 05-FEB-2024  
 Lot number: 90-24065 Exp. Date: 01-SEP-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		57.0%
Background with Soluene-350		32 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.