

### Lot specific information

**Product:** 6013181 Permafluor E+

**Mfg.Date:** 10-JUN-2024

**Batch#:** 90-24245

**Exp.Date:** 01-JAN-2027

---

#### Information

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

Test	Results
Quench parameter (tSIE)	783
Quench parameter (tSIE) with Soluene-350	567
Background with Soluene-350	45 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.**