

Lot specific information

Product: 6013231 OptiScint MP / 1L

Mfg.Date: 29-JUL-2024

Batch#: 193-24315

Exp.Date: 01-FEB-2026

Information

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

| Test | Results |
|-------------------------|---------|
| Quench parameter (tSIE) | 558 |
| Background 0-18.6 keV | 18 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.**