

## Lot specific information

Product: 6013281 OptiScint F / 1L

Mfg.Date: 10-MAR-2025

Batch#: 191-25111

Exp.Date: 01-APR-2026

## Information

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

| Test                                     | Results |
|------------------------------------------|---------|
| Quench parameter (tSIE)                  | 984     |
| Background 0-18.6 keV                    | 15 cpm  |
| Quench parameter (tSIE) with Soluene-350 | 792     |
| Background with Soluene-350              | 847 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.

revvity