

Lot specific information

Product: 6013281 OptiScint F / 1L

Mfg.Date: 11-NOV-2024

Batch#: 191-24465

Exp.Date: 01-DEC-2025

Information

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

Test	Results
Quench parameter (tSIE)	990
Background 0-18.6 keV	11 cpm
Quench parameter (tSIE) with Soluene-350	775
Background with Soluene-350	3761 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.**

