

Lot specific information

Product: 6013281 OptiScint F / 1L

Mfg.Date: 29-APR-2024

Batch#: 191-24185

Exp.Date: 01-JUN-2025

Information

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

Test	Results
Quench parameter (tSIE)	983
Background 0-18.6 keV	16 cpm
Quench parameter (tSIE) with Soluene-350	779
Background with Soluene-350	2492 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.**

