

Lot specific information

Product: 6013286 OptiScint F / 1x5L

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