

Lot specific information

Product: 6013286 OptiScint F / 1x5L

Mfg.Date: 26-MAY-2025

Batch#: 191-25221

Exp.Date: 01-JUN-2026

Information

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

Test	Results
Quench parameter (tSIE)	988
Background 0-18.6 keV	13 cpm
Quench parameter (tSIE) with Soluene-350	767
Background with Soluene-350	1547 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.