

Lot specific information

Product: 6013286 OptiScint F / 1x5L

Mfg.Date: 09-OCT-2023

Batch#: 191-23415

Exp.Date: 01-NOV-2024

Information

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

Test	Results
Quench parameter (tSIE)	997
Background 0-18.6 keV	16 cpm
Quench parameter (tSIE) with Soluene-350	794
Background with Soluene-350	1099 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.**