

Lot specific information

Product: 6013286 OptiScint F / 1x5L

Mfg.Date: 08-JUL-2024

Batch#: 191-24285

Exp.Date: 01-AUG-2025

Information

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

Test	Results
Quench parameter (tSIE)	977
Background 0-18.6 keV	18 cpm
Quench parameter (tSIE) with Soluene-350	762
Background with Soluene-350	2613 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.**