

Lot specific information

Product: 6013791 OptiScint FLOW / 1L

Mfg.Date: 15-APR-2024

Batch#: 199-24165

Exp.Date: 01-NOV-2025

Information

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

Test	Results
Quench parameter (tSIE)	503
Background 0-18.6 keV	12 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.**