

Lot specific information

Product: 6013529 Flo-Scint II

Mfg.Date: 09-SEP-2024

Batch#: 82-24375

Exp.Date: 01-APR-2026

Information

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

Test	Results
Quench parameter (tSIE)	647
Background 0-18.6 keV	16 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.**