

Lot specific information

Product: 6013167 Insta-Fluor Plus

Mfg.Date: 25-NOV-2024

Batch#: 136-24485

Exp.Date: 01-JUN-2026

Information

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

Test	Results
Quench parameter (tSIE)	853
Background 0-18.6 keV	15 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.