

Lot specific information

Product: 6013149 Filter-Count

Mfg.Date: 15-JAN-2024

Batch#: 40-24035

Exp.Date: 01-AUG-2026

Information

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

Test	Results
Quench parameter (tSIE)	665
Background 0-18.6 keV	15 cpm
Quench parameter (tSIE) with filter	616
Solubilisation speed filter	5 min

**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.**