

Lot specific information

Product: Filter-Count
 Lot number: 40-23195

Mfg. Date: 08-MAY-2023
 Exp. Date: 01-DEC-2025

Information

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

Test	Temperature	Results
Appearance		Clear blue fluorescent liquid
Quench parameter		108.0%
Background		14 cpm
Quench parameter with filter		100.0%
Water Holding Capacity	27 °C	≥ 1.5 ml clear
	5 °C	≥ 1.2 ml clear
0.15 M NaCl Holding Capacity	27 °C	≥ 2.2 ml clear
	5 °C	≥ 1.6 ml clear
Solubilisation speed filter		within 15 minutes

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.