

Lot specific information

Product: Flo-Scint II
 Lot number: 82-23365

Mfg. Date: 04-SEP-2023
 Exp. Date: 01-APR-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		100.0%
Background		14 cpm
Water Holding Capacity	27 °C	≥ 4.5 ml clear
	5 °C	≥ 1.2 ml clear
Acetonitril/water (70/30 v/v) Holding Capacity	22 °C	≥ 2.6 ml clear
1.0 M NaCl Holding Capacity	22 °C	≥ 1.5 ml clear

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.