

Lot specific information

Product: 6003038 Soluene-350

Mfg.Date: 27-MAY-2024

Batch#: 24-24225

Exp.Date: 01-DEC-2025

Information

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

Test	Results
Chlorine content	1.2 %
Normality (mg eq OH-/ g)	0.60
Quench parameter (tSIE)	570
Background 0-18.6 keV	20.00 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.**