

**Lot specific information**

**Product:** 6003038 Soluene-350

**Mfg.Date:** 28-OCT-2024

**Batch#:** 24-24445

**Exp.Date:** 01-MAY-2026

-----  
**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)	0.61
Quench parameter (tSIE)	561
Background 0-18.6 keV	16.00 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.**