

**Lot specific information**

**Product:** 6003038 Soluene-350

**Mfg.Date:** 05-AUG-2024

**Batch#:** 24-24325

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

-----  
**Information**

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

Test	Results
Chlorine content	1.1 %
Normality (mg eq OH-/ g)	0.60
Quench parameter (tSIE)	536
Background 0-18.6 keV	18.00 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.**