

### Lot specific information

**Product:** 6003005 Hydroxide of Hyamine 10-X

**Mfg.Date:** 02-OCT-2023

**Batch#:** 2-23401

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

### Information

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

Test	Results
Chlorine content	0.36 %
Normality (mg eq OH-/ g)	1.01
Quench parameter (tSIE)	347
Quench parameter Toluene scint. (tSIE)	791

**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.**