

**Lot specific information**

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

**Mfg.Date:** 10-JUN-2024

**Batch#:** 2-24245

**Exp.Date:** 01-JAN-2027

-----  
**Information**

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

Test	Results
Chlorine content	0.26 %
Normality (mg eq OH-/ g)	1.04
Quench parameter (tSIE)	344
Quench parameter Toluene scint. (tSIE)	802

**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.**

