

Lot specific information

Product: 6003005 Hydroxide of Hyamine 10-X

Mfg.Date: 04-NOV-2024

Batch#: 2-24455

Exp.Date: 01-JUN-2027

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)	1.02
Quench parameter (tSIE)	336
Quench parameter Toluene scint. (tSIE)	794

**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.**

