

**Lot specific information**

**Product:** 6013391 Insta-Gel Plus

**Mfg.Date:** 24-JUN-2024

**Batch#:** 95-24265

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

-----  
**Information**

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

Test	Results
Quench parameter (tSIE)	608
Background 0-18.6 keV	16 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.**