

Lot specific information

Product: Insta-Gel Plus
 Lot number: 95-23371

Mfg. Date: 11-SEP-2023
 Exp. Date: 01-APR-2026

Information

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

| Test | Temperature | Results |
|-------------------------|-------------|-------------------------------|
| Appearance | | Clear blue fluorescent liquid |
| Quench parameter | | 97.0% |
| Background | | 15 cpm |
| 50% water gel-formation | 5 °C | Clear rigid gel |
| Plasma Holding Capacity | 20 °C | 0.1 – 0.5 ml clear |

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.