

## Lot specific information

**Product:** 6NE9571 High Efficiency Mineral Oil Scintillator**Mfg.Date:** 10-FEB-2025**Batch#:** 137-25071**Exp.Date:** 01-SEP-2027

## Information

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

Test	Results
Quench parameter (tSIE)	685
Background 0-18.6 keV	17 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.