

Lot specific information

Product: 6013109 Monophase S

Mfg.Date: 05-FEB-2024

Batch#: 6-24065

Exp.Date: 01-SEP-2026

Information

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

Test	Results
Quench parameter (tSIE)	569
Background 0-18.6 keV	14 cpm
Background with water	19 cpm
Background Combustion	13 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.**

