

Lot specific information

Product: Monophase S
 Lot number: 6-23385

Mfg. Date: 18-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		91.0%
Background		15 cpm
Background with water		14 cpm
Background Combustion		12 cpm
Water Holding Capacity	27 °C	≥ 4.5 ml clear
	5 °C	≥ 3.2 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.