

Lot specific information

Product:	Ultima Gold MV	Mfg. Date:	05-JUN-2023
Lot number:	80-23235	Exp. Date:	01-JAN-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		104.0%
Background		16 cpm
Water Holding Capacity	27 °C	≥ 2.0 ml clear
	15 °C	≥ 1.5 ml clear
0.01 M PBS Holding Capacity	27 °C	≥ 4.5 ml clear
	15 °C	≥ 4.0 ml clear
0.15 M NaCl Holding Capacity	27 °C	≥ 4.0 ml clear
	15 °C	≥ 3.5 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.