

Lot specific information

Product: Ultima Gold XR

Mfg. Date: 09-OCT-2023

Lot number: 79-23415

Exp. Date: 01-MAY-2025

Information

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

Test	Temperature	Results
Appearance		Clear blue fluorescent liquid
Quench parameter		88.0%
Background		16 cpm
Water Holding Capacity	27 °C	2 - 10 ml clear
	15 °C	2 - 10 ml clear
0.01 M PBS Holding Capacity	27 °C	2 - 10 ml clear
	15 °C	2 - 10 ml clear
0.5 M (NH ₄) ₂ SO ₄ Holding Capacity	27 °C	≥ 1.1 ml clear
	5 °C	≥ 1.3 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.**