



# DISCOVERY

L I F E S C I E N C E S

## CERTIFICATE OF ANALYSIS

**Product Name:** Universal Cryopreservation Recovery Medium (UCRM™)  
50 mL

**Catalog Number:** 4.81015      **Storage Conditions:** 4°C  
**Lot Number:** 2408291      **Date Released:** September 11, 2024  
**Product Brand:** Gentest®      **Expiration Date:** May 16, 2026

ATTRIBUTES	SPECIFICATION	RESULTS
Appearance	Translucent pink liquid in 50mL tube	CONFORMS
pH	7.4 ± 0.2	PASS
7-day Sterility	Sterile	PASS

Quality Assurance

12 September 2024

Date