



DISCOVERY

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

CERTIFICATE OF ANALYSIS

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

Catalog Number: 4.81074 **Storage Conditions:** 4°C
Lot Number: 2402272 **Date Released:** March 05, 2024
Product Brand: Gentest® **Expiration Date:** January 08, 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

18 March 2024

Date