



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: 2408160 **Date Released:** September 3, 2024
Product Brand: Gentest® **Expiration Date:** August 27, 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

10 September 2024

Date