



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: 2408290 **Date Released:** September 11, 2024
Product Brand: Gentest® **Expiration Date:** July 2, 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