



DISCOVERY

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

CERTIFICATE OF ANALYSIS

Product Name: Universal Primary Cell Plating Medium (UPCM™) 50 mL

Catalog Number: 4.81016 **Storage Conditions:** 4°C
Lot Number: 2403270 **Date Released:** April 8, 2024
Product Brand: Gentest® **Expiration Date:** April 1, 2026

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

Quality Assurance

15 March, 2024

Date