



DISCOVERY

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

CERTIFICATE OF ANALYSIS

Product Name: HIM™: Hepatocyte Induction Medium 50mL (with Supplements)

Catalog Number: 4.81018 **Storage Conditions:** 4°C
Lot Number: 2403273 **Date Released:** April 18, 2024
Product Brand: Gentest® **Expiration Date:** July 11, 2024

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

Quality Assurance

03 May 2024

Date