



DISCOVERY

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

CERTIFICATE OF ANALYSIS

Product Name: RHIM™: Rodent Hepatocyte Induction Medium, 50 mL (with Supplements)

Catalog Number: 4.81037 **Storage Conditions:** 4°C
Lot Number: 2404154 **Date Released:** April 23, 2024
Product Brand: Gentest® **Expiration Date:** July 16, 2024

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

Quality Assurance

09 May 2024

Date