



# 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:** 2409204      **Date Released:** October 16, 2024  
**Product Brand:** Gentest®      **Expiration Date:** January 9, 2025

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

Quality Assurance

October 21 2024

Date