


## CERTIFICATE OF ANALYSIS

Product Name: **CHPM™ Cryopreserved Hepatocyte Plating Medium**  
 Product Number: 4.70002 (CM9000)  
 Product Brand: Gentest®  
 Storage Temperature: 4-8°C DO NOT FREEZE.

Lot Number:	Date of Manufacture:	Expiration Date (2 years from date of manufacture):	Initials and Date of CoA Preparer:
2408081	August 12, 2024	August 12, 2026	KJG 8/19/24

TEST	SPECIFICATION	LOT RESULTS	Initials and Date of Inspector
Appearance – translucent liquid in 50 ml tube	Pass/fail	Pass	KJG 8/19/24
PH	7.40 ± 0.2	7.39	KJG 8/19/24
Sterility Test: tested in medium for 7 days in incubator to indicate no visual contamination	Pass/Fail	Pass	KJG 8/19/24
Inclusions	None	None	KJG 8/19/24
Leakage	None	None	KJG 8/19/24
Cap and Tube	Intact	Intact	KJG 8/19/24

**Quality Review and Release:**

Name (Printed) and title:	Edward Francis Senior Quality Specialist
Signature:	
Date:	23 August 2024