


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:
2410112	10/22/24	10/22/26	KJG 11/4/24

TEST	SPECIFICATION	LOT RESULTS	Initials and Date of Inspector
Appearance – translucent liquid in 60 ml PETG bottle	Pass/fail	Pass	KJG 11/4/24
PH	7.40 ± 0.2	7.40	KJG 11/4/24
Sterility Test: tested in medium for 7 days in incubator to indicate no visual contamination	Pass/Fail	Pass	KJG 11/4/24
Inclusions	None	None	KJG 11/4/24
Leakage	None	None	KJG 11/4/24
Cap and Tube	Intact	Intact	KJG 11/4/24

Quality Review and Release:

Name (Printed) and title:	Edward Francis, Sr. Quality Specialist
Signature:	
Date:	November 6, 2024