



## CERTIFICATE OF ANALYSIS

Product Name: **CellAffix™ - Collagen Coated Tissue Culture Plate (96-Well Plate)**

Product Number: 4.71021 (CM1096)

Product Brand: Gentest®

Storage Temperature: 4-8°C

TEST	SPECIFICATION	LOT # 2407174 RESULTS
Appearance	clear multi-well tissue culture plate	Conforms
Coating	Collagen Coating	Conforms
Shelf Life	5 years	August 26, 2029
QC acceptance date	7 days clearance	September 4, 2024

Signature:

A handwritten signature in black ink, appearing to read "Edward Francis".

Printed Name:

Edward Francis

Title:

Sr. Quality Specialist

Date:

05 September 2024