



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 # 2407175 RESULTS
Appearance	clear multi-well tissue culture plate	Conforms
Coating	Collagen Coating	Conforms
Shelf Life	5 years	August 28, 2029
QC acceptance date	7 days clearance	September 6, 2024

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

Signature:

Printed Name: Edward Francis

Title: Sr. Quality Specialist

Date: 10 September 2024