



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 # 2406070 RESULTS
Appearance	clear multi-well tissue culture plate	Conforms
Coating	Collagen Coating	Conforms
Shelf Life	5 years	June 10, 2029
QC acceptance date	7 days clearance	June 18, 2024

Signature:

A handwritten signature in black ink, appearing to read "E. Francis", written over a light blue background.

Printed Name:

Edward Francis

Title:

Sr. Quality Specialist

Date:

21 June 2024