



DISCOVERY

L I F E S C I E N C E S

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 # 2407171 RESULTS
Appearance	clear multi-well tissue culture plate	Conforms
Coating	Collagen Coating	Conforms
Shelf Life	5 years	August 19, 2029
QC acceptance date	7 days clearance	August 28, 2024

Signature:

Alexa Silies

Printed Name:

Alexa Silies

Title:

Quality Assurance Specialist

Date:

08/30/2024