



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 5 pack)

Catalog Number: 4.71008 **Storage Conditions:** 4°C
Lot Number: 2408294 **Date Released:** September 12, 2024
Product Brand: Gentest® **Expiration Date:** September 03, 2029

ATTRIBUTES	SPECIFICATION	RESULTS
Appearance	Clear multi-well tissue culture plate	CONFORMS
Coating	Collagen Coating	CONFORMS
7-day Sterility	PASS	PASS

Quality Assurance

18 September, 2024

Date