



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: 2403182 **Date Released:** April 5, 2024
Product Brand: Gentest® **Expiration Date:** March 28, 2029

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

Quality Assurance

22 April 2024

Date