



**G-Biosciences** ♦

♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

## Certificate of Analysis / Data Sheet

**Product:** Well-Coat Protein A

**Catalog No:** 786-770

**Release Date:** 11/27/19

**Size:** 5 plates

**Retest Date:** 11/27/25

**Lot No:** 192711

**Storage:** 4°C

### Specifications

**Appearance:** Black 96-well plates

Test	Specification	Result
Visual	Plates clear, colorless, free of foreign matter, and scratches	PASS
ELISA(Average CV)	< 10%	PASS
ELISA(signal to noise ratio)	> 7.0	PASS
ELISA(plate to plate CV)	≤ 15%	PASS

This document has been produced electronically and is valid without signature

**Date of Print:** December 4, 2019

**G-Biosciences/Geno Technology, Inc.**

Rev 10/25/18 BP



**think proteins! think G-Biosciences** [www.GBiosciences.com](http://www.GBiosciences.com)