



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Well-Coat Protein A

Catalog No: 786-731

Release Date: 11/27/19

Size: 5 plates

Retest Date: 11/27/25

Lot No: 192711

Storage: 4°C

Specifications

Appearance: Clear strip 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