



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: C18 Spin Columns

Catalog No: 786-931

Release Date: 08/16/2021

Size: 50 columns

Retest Date: 08/16/2026

Lot No: 211608

Storage: Room Temp.

Specifications

CAS Number: N/A

Appearance: White, dry spherical beads

Test	Specification	Result
Particle Size	15µm	PASS
Pore Size	300Å	PASS
Quantity per column	10mg	PASS

This document has been produced electronically and is valid without signature

Date of Print: December 21, 2021

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



think proteins! think G-Biosciences www.GBiosciences.com