



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: SpinOUT™ GT-1200

Catalog No: 786-173

Release Date: 08/11/2021

Size: 10 x 3ml columns

Retest Date: 08/11/2026

Lot No: 211108

Storage: 4°C in 20% ethanol

Specifications

CAS Number: N/A

Appearance: White resin in clear solution

Test	Specification	Result
Functional	Performs as per corresponding and/or companion protocol	PASS

This document has been produced electronically and is valid without signature

Date of Print: January 20, 2022

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



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