



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: SpinOUT™ GT-600, 1ml

Catalog No: 786-720

Release Date: 03/08/2022

Size: 5 x 10/bag

Retest Date: 03/08/2027

Lot No: 220803

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: May 11, 2022

G-Biosciences/Geno Technology, Inc.

Rev 01/20/22 DB



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