



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Concanavalin A (Con A) Agarose

Catalog No: 786-217

Release Date: 06/25/2020

Size: 25 ml resin

Retest Date: 06/25/2023

Lot No: 202506

Storage: 4°C

Specifications

CAS Number: N/A

Appearance: 50% slurry. White agarose beads in clear solution.

Test	Specification	Result
Ligand Density	10-16mg Con A/ml resin	PASS
Capacity	20-50mg thyroglobulin/ml resin	PASS
Bead Structure	6% agarose	PASS
pH Stability	4.0-9.0	PASS

This document has been produced electronically and is valid without signature

Date of Print: January 28, 2022

G-Biosciences/Geno Technology, Inc.

Rev 01/24/22 DB



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