



G-Biosciences ♦

♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Copper Chelating Resin

Catalog No: 786-285

Release Date: 05/17/2023

Size: 10 ml

Retest Date: 05/07/2029

Lot No: 230905

Storage: 4 °C

Specifications

Appearance: Blue resin in clear solution
Bead Structure: 6% cross-linked agarose

Test	Specification	Result
Binding capacity	> 50 mg his tagged protein/ml resin	PASS
Ligand density	20 – 40 μ moles Cu^{2+} /ml resin	PASS

This document has been produced electronically and is valid without signature

Date of Print: June 9, 2023

G-Biosciences/Geno Technology, Inc.

Rev 03/28/2023 DB



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