



Certificate of Analysis / Data Sheet

Product: Ni-NTA Resin

Catalog No: 786-939 to 786-945, & 786-1547

Lot No: 191710

Size: 10mL, 100mL, 500mL, 2 x
500mL, 25mL, 25 x 0.2mL, 5 x
1mL, 5 x 3mL

Release Date: 10/17/19

Retest Date: 10/17/24

Storage: 4°C in 20% ethanol. DO NOT FREEZE.

Specifications

Matrix: Cross-linked agarose, 6%
Particle Size: Spherical, diameter 50-160µm
Max Pressure (MPa): 0.02
pH Stability: 3-12

Test	Specification	Result
Ni load/mL resin	≥16µmol	PASS
Visual inspection	Green Slurry	PASS
Slurry percentage	50% resin volume	PASS

This document has been produced electronically and is valid without signature

Date of Print: November 18, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

