



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

Certificate of Analysis

Product: Ni-NTA Resin
Catalog #: 786-939, 786-940, 786-941, 786-942, 786-943, 786-944, 786-945
Lot#: 183004
Matrix: Cross-linked agarose, 6%
Particle Size: Spherical, diameter 50-160µm
Max Pressure (MPa): 0.02
pH Stability: 3-12
Storage Condition: 4-30°C, 20% ethanol

Test(s)	Specifications	Result
Ni Load /ml Resin	≥16 µmol	Passed
Visual Inspection	Green slurry	Passed
Slurry Percentage	50% resin volume	Passed

Storage: Store at 4°C. DO NOT FREEZE.

This document has been produced electronically and is valid without signature

Date of Print: May 17, 2018

For Research Use Only

G-Biosciences / Genotech
9800 Page Avenue
St. Louis, MO 63132, USA
Ph: 314-991-6034, Fax: 314-991-1504



think proteins! think G-Biosciences!

www.GBiosciences.com