



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

Certificate of Analysis / Data Sheet

Product: Alum Adjuvant

Catalog No: 786-1215

Size: 50 ml

Lot No: 231505

Storage: 4 °C

Specifications

Test Procedure

White colloidal suspension	Verified	Yes / No
Aluminum content (~10mg/ml)	Verified	Yes / No
pH (6.5-7.0)	Verified	Yes / No
Sterility	Verified	Yes / No

This document has been produced electronically and is valid without signature

Date of Print: June 29, 2023

G-Biosciences/Geno Technology, Inc.

9800 Page Avenue, St. Louis, MO 63132, USA

Ph: 314-991-6034, Fax: 314-991-1504

Rev 02/13/17 CCH



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