



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

## **CERTIFICATE OF ANALYSIS**

**Product Name:** DNA *OUT*<sup>™</sup>

**Catalog #** 786-74

**Lot #:** 210812

**Size:** 250ml Spray Bottle

### **Tests**

DNA <i>OUT</i> appearance	Clear Liquid	Passed	Yes / <del>No</del>
---------------------------	--------------	--------	---------------------

pH Test		Passed	Yes / <del>No</del>
---------	--	--------	---------------------

**Storage Temp:** Store at Room Temp

*This document has been produced electronically and is valid without a signature*

*For Research Use Only*

**Print Date:** December 27, 2021

DNAOut C of A/Common#3  
Rev 02.19.13/IA



think proteins! think G-Biosciences!

[www.GBiosciences.com](http://www.GBiosciences.com)