



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

Certificate of Analysis

Product: Alamar Blue Cell Viability Reagent

Catalog #: 786-923

Lot #: 190212

Re-test date: 12-2024

Tests:

Appearance	Dark blue solution with no visible ppt.
Release Method	Absorbance 0.74-0.83 AU
Microbiology testing	No growth at 34°C to 36°C for minimum of 72 hours

Storage: Store at 4°C *for Research Use Only*

This document has been produced electronically and is valid without signature

Date of Print: December 19, 2019

G-Biosciences / Genotech

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

Alamar blue/Com#3
Rev 04.14.17CCH



think proteins! think G-Biosciences!

www.GBiosciences.com