



## Certificate of Analysis

**Product:** Phenol Red Solution 0.5%

**Catalog #:** 786-563

**Size:** 250ml

**Lot#:** 222401

**Re-test date:** 01/2027

### **Test Specification:**

Appearance- Dark Red Color	Passed	Yes / <del>No</del>
pH Test at RT (7.0 – 7.6)	Passed	Yes / <del>No</del>
Osmolality with NaHCO <sub>3</sub> (291 -321 MOSM/kg)	Passed	Yes / <del>No</del>

**Storage Temp:** Store at Room Temp.

This document has been produced electronically and is valid without signature

**Date of Print:** February 4, 2022

### **G-Biosciences / Genotech**

9800 Page Avenue

St. Louis, MO 63132, USA

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

Phenol Red Solution 0.5% C of A/Com#3  
Rev 08.24.12/MM

