



Certificate of Analysis

Product: Phenol Red Solution 0.5%

Catalog #: 786-563

Size: 250ml

Lot#: 203006

Re-test date: 06/2025

Test Specification:

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

Storage Temp: Store at Room Temp.

This document has been produced electronically and is valid without signature

Date of Print: August 5, 2020

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

