



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:** CHAPS 5% Solution (Filtered)

**Catalog #:** DG099

**Size:** 500ml

**Lot#:** 161804

**Appearance:** Clear liquid

**Tests:**

|                  |                       |        |
|------------------|-----------------------|--------|
| Conductivity     | 2.2ppm / 4.38 $\mu$ S | PASSED |
| pH               | 5.20                  | PASSED |
| Specific gravity | 1.01                  | PASSED |

**Shelf Life:** May 31, 2017

**Storage:** Store at 4<sup>0</sup>C.

This document has been produced electronically and is valid without signature

*For Research Use Only*

**Date of Print:** April 26, 2016

**G-Biosciences / Genotech**

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

CHAPS, 5% Solution (Filtered) C of A/Com3  
Rev 09.17.15/IA



think proteins! think G-Biosciences!

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