



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

Certificate of Analysis

Product: CHAPS 5% Solution (Filtered)

Catalog #: DG099

Size: 500ml

Lot#: 162106

Appearance: Clear liquid

Tests:

Conductivity	2.2ppm / 4.38 μ S	PASSED
pH	5.20	PASSED
Specific gravity	1.01	PASSED

Shelf Life: May 31, 2017

Storage: Store at 4⁰C.

This document has been produced electronically and is valid without signature

For Research Use Only

Date of Print: September 9, 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