



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 #: DG097

Size: 50ml

Lot#: 101011

Appearance: Clear liquid

Tests:

Conductivity	3.07ppm / 6.12 μ S	PASSED
pH	4.6	PASSED
Specific gravity	1.04	PASSED

Storage: Store at 4⁰C.

For Research Use Only

This document has been produced electronically and is valid without signature

Date of Print: December 2, 2010

G-Biosciences / Genotech

92 Weldon Parkway, St. Louis, MO 63043, USA
Fax: 314-991-1504

CHAPS 5% Solution (filtered) C of A/Com3
Rev 11.19.10/IA



think proteins! think G-Biosciences!

www.GBiosciences.com