



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; [3-(3-Cholamidopropyl) dimethylammonio-1-propanesulfonate]  
**Catalog #:** DG049  
**Size:** 1g  
**Lot#:** 102707  
**Appearance:** White Powder  
**Mol. Formula:**  $C_{32}H_{58}N_2SO_7$   
**Mol Weight:** 614.89  
**CAS#:** [75621-03-3]

### Tests:

Assay (from N)	98.9 %	PASSED
Solubility (20%, Water)	Colorless	PASSED
Nitrogen (Kjeldahl)	4.5%	PASSED
Moisture	0.89%	PASSED
TLC One Spot (Silica gel: methanol/NH <sub>4</sub> OH = 95/5)		PASSED
Conductivity (0.1M in H <sub>2</sub> O)	4.14µs/cm	PASSED

**Storage:** Store at Room Temp. Protect from Moisture.

This document has been produced electronically and is valid without signature

**Date of Print:** April 6, 2011

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

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

CHAPS DG050 25g C of A/Com3  
Rev 12.18.08-SA/MM/IA



think proteins! think G-Biosciences!

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