



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:** CHAPSO; [3-(3-Cholamidopropyl) dimethylammonio]-2-hydroxy-1-propanes

**Catalog #:** DG052

**Size:** 1g

**Lot#:** 091211

**Appearance:** White Powder

**Mol Weight:** 630.9 daltons

**CAS#:** [82473-24-3]

### Tests:

Infrared Spectrum	Matches reference spectrum	PASSED
Solubility (10%, Water)	Colorless	PASSED
TLC One Spot (Silica gel: methanol/NH <sub>4</sub> OH = 95/5)		PASSED
Conductivity (0.5M in H <sub>2</sub> O)	12 micromhos	PASSED

**Storage:** Store at 4° C. Protect from Moisture.

This document has been produced electronically and is valid without signature

**Date of Print:** November 12, 2009

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

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

CHAPSO DG052 1g C of A/Com3  
Rev 12.18.08-SA/MM/IA



**think proteins! think G-Biosciences!**

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