



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

Certificate of Analysis

Product: CHAPS; [3-(3-Cholamidopropyl) dimethylammonio-1-propanesulfonate]
Catalog #: DG050
Size: 5g
Lot#: 092008
Appearance: White Powder
Mol. Formula: C₃₂H₅₈N₂SO₇
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 ₄ OH = 95/5)		PASSED
Conductivity (0.1M in H ₂ 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: August 24, 2009

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