



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 #: DG051
Size: 25g
Lot#: 101407
Appearance: White Powder
Mol. Formula: C₃₂H₅₈N₂SO₇
Mol Weight: 614.89
CAS#: [75621-03-3]

Tests:

Assay (from N)	99.91 %	PASSED
Solubility (10% in Water, 20°C)	Colorless	PASSED
Nitrogen (Kjeldahl)	4.55%	PASSED
Moisture	0.71%	PASSED
TLC One Spot (Silica gel: methanol/NH ₄ OH = 95/5)		PASSED
Conductivity (0.1M in H ₂ O)	6.15µs/cm	PASSED

Storage: Store at Room Temp. Protect from Moisture.

This document has been produced electronically and is valid without signature

For Research Use Only

Date of Print: December 3, 2010

Genotech / G-Biosciences

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

CHAPS DG051 25g C of A/Com3/Rev 12.18.0-SA/MM/IA



think proteins! think G-Biosciences!

www.GBiosciences.com