



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

Certificate of Analysis

Product: CHAPSO; [3-(3-Cholamidopropyl) dimethylammonio-1-propanesulfonate]

Catalog #: DG053

Size: 5g

Lot#: 120202

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 ₄ OH = 95/5)	PASSED
Conductivity (0.5M in H ₂ 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: February 2, 2012

G-Biosciences / Genotech

92 Weldon Parkway, St. Louis, MO 63043, USA

Fax: 314-991-1504

CHAPSO DG053 5g C of A/12.03.09-IA/Common3



think proteins! think G-Biosciences!

www.GBiosciences.com