



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]-2-hydroxy-1-propanesulfonate

Catalog #: DG053

Size: 5g

Lot#: 121207

Appearance: White powder

Mol Weight: 630.89 Daltons

CAS#: [82473-24-3]

Tests:

Infrared Spectrum - Matches reference spectrum	PASSED
Iron (Fe) <1ppm	PASSED
Conductivity <43 µmho/cm	PASSED

Storage: Store at 4° C. Protect from Moisture and Light.

This document has been produced electronically and is valid without signature

Date of Print: July 13, 2012

G-Biosciences / Genotech

9800 Page Avenue, St. Louis, MO 63132, USA

Fax: 314-991-1504

CHAPSO DG053 5g C of A/07.12.12/MM



think proteins! think G-Biosciences!

www.GBiosciences.com