



### Certificate of Analysis

**Product:** CHAPS; [3-(3-Cholamidopropyl) dimethylammonio-1-propanesulfonate]  
**Catalog #:** DG050  
**Size:** 5g  
**Lot#:** 203011  
**Appearance:** White Powder  
**Mol. Formula:** C<sub>32</sub>H<sub>58</sub>N<sub>2</sub>SO<sub>7</sub>  
**Mol Weight:** 614.89  
**CAS#:** [75621-03-3]  
**Re-test date:** 11/2026

**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 <sub>4</sub> OH = 95/5)		PASSED
Conductivity (0.1M in H <sub>2</sub> 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 21, 2020

**Genotech / G-Biosciences**

9800 Page Avenue  
St. Louis, MO 63132, USA  
Fax: 314-991-1504

CHAPS DG051 25g C of A/Com3 Rev 05.22.12/MM

