



Certificate of Analysis / Data Sheet

Product: CHAPSO

Catalog No: DG052

Release Date: 2/4/19

Size: 1g

Retest Date: 1/17/23

Lot No: 191701

Storage: 4°C

Specifications

CAS Number: 82473-24-3

Appearance: White powder

Test	Specification	Result
Solubility	10% clear aqueous solution	Meets Specification
TLC	One spot	Meets Specification
Conductivity	0.5M <100micromhos	12micromhos
Absorbance (10% water)	A400 <0.010% A340 <0.020%	Meets Specification
Karl Fischer	<6%	2.8%
Infrared Spectrum	Matches reference spectrum	Meets Specification

This document has been produced electronically and is valid without signature

Date of Print: December 20, 2019

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP

