



Certificate of Analysis / Data Sheet

Product: CMC-535™ Detergent Assay

Catalog No: DG535

Lot No: 201106

Components: Fluorescent Dye [100X]

Release Date: 6/11/2020

Reagent 1

Retest Date: 7/11/2021

Reagent 2 [100X]

Storage: 4°C

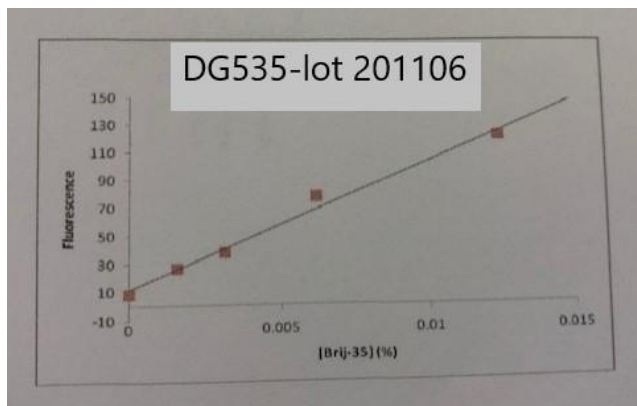
Specifications

Appearance: Fluorescent Dye [100X] yellow clear solution

Appearance: Reagent 1 clear solution

Appearance: Reagent 2 [100X] colorless clear solution

Test Result



This document has been produced electronically and is valid without signature

Date of Print: January 28, 2022

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP



