



## Certificate of Analysis / Data Sheet

**Product:** SDS (Sodium dodecyl sulfate)

**Catalog No:** DG093

**Release Date:** 9/20/2021

**Size:** 500g

**Retest Date:** 10/30/2023

**Lot No:** 211509

**Storage:** Room temp.

### Specifications

**CAS Number:** 151-21-3

**Appearance:** Fine white to off-white powder

Test	Specification	Result
Abs.@260nm (3%, Water)	$\leq 0.1$	0.0162
Abs.@280nm (3%, Water)	$\leq 0.1$	0.0093
Chloride	$\leq 0.1\%$	0.01%
Copper	$\leq 0.0005\%$	$< 0.0005\%$
DNase	NONE / Fail	NONE
Heavy Metals (as Pb)	$\leq 0.0005\%$	$< 0.0002\%$
Identification	PASS / Fail	PASS
Iron	$\leq 0.0001\%$	$< 0.0001\%$
Phosphate	$\leq 0.0005\%$	$< 0.0001\%$
Protease	NONE / Fail	NONE
Purity	$> 99.0\%$	99.8%
RNase	NONE / Fail	NONE
Solubility (10%, Water)	PASS / Fail	PASS

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

