



Certificate of Analysis / Data Sheet

Product: L-Ascorbic Acid**Catalog No: RC-022****Release Date: 10/9/20****Size: 100g****Retest Date: 10/9/23****Lot No: 200910****Storage: Room Temp.****Specifications****CAS Number: 50-81-7****Appearance: White crystalline powder**

Test	Specification	Result
Assay	99.0-100.5%	99.78%
Identification	Positive Reaction	PASS
Melting Point	About 190°C	190.9°C
pH (2% solution)	2.4-2.8	2.43
pH (5% solution)	2.1-2.6	2.36
Solubility in water 80%	0.33g/ml	PASS
Clarity of Solution	Clear	PASS
Color of Solution	BY ₇ Maximum	PASS
Loss on Drying	≤ 0.4%	< 0.4%
Residual Solvents	Meets specs.	PASS
Specific Optical Rotation	+20.5° to +21.5°	+21.08°
Sulfate Ash	≤ 0.1%	0.1%
Copper	≤ 5ppm	< 5ppm
Heavy Metals	≤ 10ppm	< 5ppm
Iron	≤ 2ppm	< 2ppm
Lead	≤ 2mg/kg	< 2mg/kg
Cadmium	≤ 1mg/kg	< 1mg/kg



Foreign Matter	$\leq 2.0\text{mm}$	$< 2.0\text{mm}$
Oxalic Acid	$\leq 0.2\%$	$< 0.2\%$
Arsenic	$\leq 3\text{ppm}$	$< 3\text{ppm}$
Mercury	$\leq 0.1\text{mg/kg}$	$< 0.1\text{mg/kg}$
Zinc	$\leq 0.25\text{ppn}$	$< 0.25\text{ppm}$
Salmonella	Negative	PASS
E. Coli	Negative	PASS
Pseudomonas	Negative	PASS
Staphylococcus	Negative	PASS
Total Plate Count	$\leq 100\text{cfu/g}$	$< 100\text{cfu/g}$
Yeast and Mold	$\leq 100\text{cfu/g}$	$< 100\text{cfu/g}$
BSE/TSE	BSE/TSE Free	BSE/TSE Free

<p>This document has been produced electronically and is valid without signature</p>
--

Date of Print: October 20, 2020

G-Biosciences/Geno Technology, Inc.

Rev 10/25/18 BP