



G-Biosciences, St Louis, MO. USA ♦ 1-800-628-7730 ♦ 1-314- 991-6034 ♦ technical@GBiosciences.com

Certificate of Analysis

Product: L-Lysine monohydrate

Catalog #: RC-066

Size: 100g

Lot#: 122706

Appearance: White powder

Mol Weight: 164.21

Tests	Min	Max	Actual	Units	Disposition
Protease		None	None	P/F	Passed
Heavy Metals		10	<=5	ppm	Passed
TLC, Other Amino acids		0.5	<=0.5	%	Passed
Chloride		300	<=300	ppm	Passed
Sulfate		200	<=200	ppm	Passed
Residue on Ignition		0.1	0.00	%	Passed
Water	10	11.5	11.29	%	Passed
Specific Rotation	+25.5	+27.0	+26.211	degrees	Passed
pH (2.5%, Water) at 25C	9.0	10.5	10.0		Passed
Transmittance	95.0		99.60	%	Passed

Storage: Store at 4°C. Keep dry. Warm it to room temp before opening.

This document has been produced electronically and is valid without signature

Date of Print: June 27, 2012

For Research Use Only

G-Biosciences / Genotech

9800 Page Avenue
St. Louis, MO 63132, USA
Ph: 314-991-6034, Fax: 314-991-1504

Lysine C of A/Rev 6.27.12/IA



think proteins! think G-Biosciences!

www.GBiosciences.com