



655 Research Parkway
Suite 525
Oklahoma City, OK 73104
Phone: 405-271-1544
Fax: 405-271-1545
E-mail: info@hyalose.com
www.hyalose.com

Certificate of Analysis

nanoHA-5 Low Endotoxin

Cat. Number: HYA-NANO5-1

Lot No.: H-N5-070511

Appearance

Ranging from white crystalline to clear glassy solids

Purity

May contain trace amounts of other nanoHA or oligoHA sizes

Quantity

1 mg carbohydrate (as determined
by carbazole assay)

Maximal Protein Level

Not determined

Maximal Endotoxin Level

≤ 0.04 EU/mg oligosaccharide

QC Acceptance Date

May 2007

A handwritten signature in blue ink, appearing to read "M. Haller", positioned above a horizontal line.

Michael Haller, Ph.D.

A handwritten signature in blue ink, appearing to read "Wei Jing", positioned above a horizontal line.

Wei Jing, Ph.D., M.D.