



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

Select-HA™ B500

Catalogue No. HYA-B500-200

Lot No. B500-200603

Molecular Mass: 520 kDa

Appearance

Ranging from white crystalline to clear glassy solids

Quantity

200 pmole (~104 µg)

End-label

Biotin

Protein content

<0.1%

Molecular weights

Values as determined by Multi-angle Laser Light Scattering-Size Exclusion Chromatography (SEC-MALLS) are typically within ± 5%.

M_n, 518.8 kDa

M_w, 519.8 kDa

M_w/M_n, 1.002

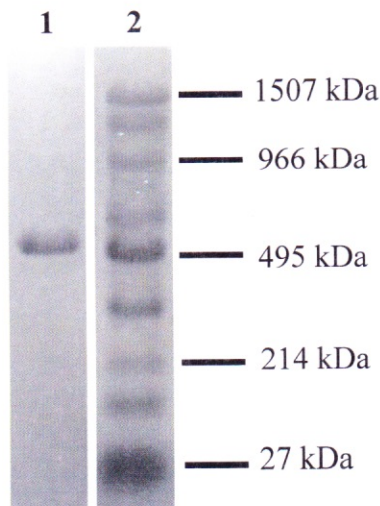


Figure 1. Electrophoretic analysis of Select-HA B500-200603.

Select-HAs were run on 0.7% agarose gel and stained with Stains-all (0.005 – 0.01% of Stains-all in 50% ethanol; overnight with gentle shaking).

lane 1, B500-200603; **lane 2**, Select-HA™ Ladder (molecular masses are indicated).

QC Acceptance Date

March 2006



Wei Jing, MD., Ph.D.



Michael Haller, Ph.D.