



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™ B250

Catalogue No. HYA-B250-200

Lot No. B250-200603

Molecular Mass: 280 kDa

Appearance

White powder

Quantity

200 pmole (~56 µg)

End-label

Biotin

Protein content

<0.1%

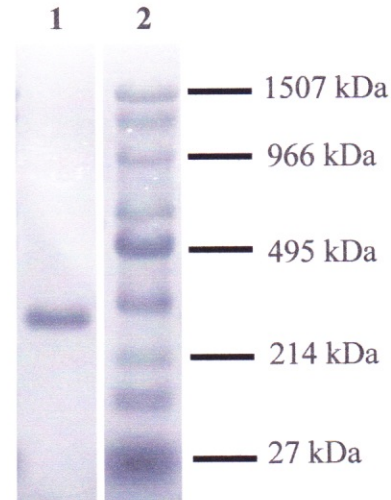


Figure 1. Electrophoretic analysis of Select-HA B250-200603.

Select-HAs were run on 0.7% agarose gel and stained with Stains-all. **lane 1**, B250-200603; **lane 2**, Select-HA™ Ladder (molecular masses are indicated).

Molecular weights

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

M_n, 279.2 kDa

M_w, 279.6 kDa

M_w/M_n, 1.002

QC Acceptance Date

March 2006


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


Michael Haller, Ph.D.