



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™ HiLadder
Catalogue No. HYA-HILAD-20
Lot No. HL201504

Appearance
Clear solid.

Quantity
~25 µg of HA.

Solubility
Soluble in water.

1648 kDa —
1138 kDa —
940 kDa —
667 kDa —
509 kDa —



Figure 1.
Electrophoretic Separation of Select-HA™ HiLadder Standards.
Five µl of reconstituted HiLadder sample was separated on a 0.7% (w/v) agarose gel by electrophoresis and stained with Stains-all.

Number-average molecular mass (<i>M_n</i> , Da)	Weight-average molecular mass (<i>M_w</i> , Da)	Polydispersity (<i>M_w/M_n</i>)
1622 k	1648 k	1.016
1114 k	1138k	1.022
864 k	940 k	1.088
660 k	667 k	1.011
507 k	509 k	1.003

Molecular Masses

Values as determined by Multi-angle Laser Light Scattering-Size Exclusion Chromatography (MALLS-SEC) are typically within ± 5%. Results in the table are the average of duplicate analysis.

QC Acceptance Date
March 2015


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


Jonathan Roberts, BS