

# Echelon Biosciences Inc.

## Select-HA™

Product Name	Catalog Number	Molecular Mass
Select-HA™ 25k	HYA-0025	18-33 kDa
Select-HA™ 50k	HYA-0050	38-63 kDa
Select-HA™ 75k	HYA-0075	68-83 kDa
Select-HA™ 200k	HYA-0200	180-220 kDa
Select-HA™ 400k	HYA-0400	360-440 kDa
Select-HA™ 500k	HYA-0500	450-550 kDa
Select-HA™ 600k	HYA-0600	570-630 kDa
Select-HA™ 900k	HYA-0900	855-945 kDa
Select-HA™ 1000k	HYA-1000	950-1050 kDa
Select-HA™ 1250k	HYA-1250	1125-1375 kDa

Support: echelon@echelon-inc.com

### Description

Select-HA™ is a hyaluronic acid (HA) preparation of uniform and narrow size distribution prepared by in vitro synthesis using recombinant *Pasteurella multocida*

hyaluronan synthase<sup>1</sup>. Select-HA™ is a trademark of Hyalose LLC.

### Properties

**Size** - 1 mg. Determined by Carbazole Assay based on free acid form. Due to the existence of sodium, the actual dry weight is more than 1 mg.  
**Form** - lyophilized hyaluronan polymers as sodium salts

**Storage** - -20 °C or below. Avoid frequent freeze-thaw, aliquoting is recommended. Avoid contamination with microbes or HA-degrading enzymes.

**Reconstitution** - Carefully open vial and add desired amount of sterile water or sterile desired buffer. Ensure HA powder at the bottom or on the sides of the vial are dissolved. Allow two hours at 4 °C with periodic mixing for full rehydration. Centrifuge vial for a few seconds to collect the HA solution at the bottom of the tube.

**Molecular Mass** - see table

**Endotoxin Level** - < 0.1 EU/mg

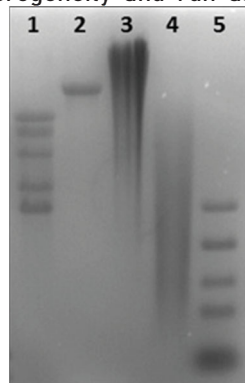
\*Please see Certificate of Analysis for lot specific information

### Background

Hyaluronic acid (HA) is a high molecular weight anionic polysaccharide (1,000-10,000 kD) composed of repeating disaccharides and is one of several glycosaminoglycan components of the extracellular matrix of connective tissue. Free HA is taken up by the liver where it is degraded and recycled. Data indicates a relationship between HA levels, local inflammation, and severity of many diseases such as hepatitis B or C, rheumatoid arthritis, liver fibrosis, etc.

### Data: Agarose Gel

Select-HA™ (lane 2) and Select-HA Ladder™ (lane 1 & 5) shows tight bands in agarose gel while other commercial HA (lane 3 & 4) have much greater size heterogeneity and run as a smear.



### References

1. Jing, W.; DeAngelis, P. L. (2004) Synchronized chemoenzymatic synthesis of monodisperse hyaluronan polymers. *J Biol Chem*, 279 (40), 42345-9.

### Related Products

Product	Catalog Number
<b>Compounds</b>	
BODIPY-HA	H-025F, H-250F, H-700F
Texas Red-HA	H-025R, H-250R, H-700R
Biotinylated Select-HA™	HYA-B50-200, HYA-B250-200, HYA-B500-200, HYA-B1000-200
Select-HA Ladder™	HYA-HILAD-20, HYA-LOLAD-20, HYA-MGLAD
nanoHA™	HYA-NAN05-1
HAase Inhibitor	B-0601
<b>HA Binding Proteins</b>	
Versican G1 Domain	G-HA01, G-HA02
<b>Assays</b>	
HA Quantification ELISAs	K-1200, K-4800
Hyaluronidase Activity ELISA	K-6000

Technical Data Sheet Rev. 4 10-12-23 - For research use only. Not intended or approved for diagnostic or therapeutic use.



Echelon Biosciences Inc.

Ph: 866-588-0455

Fax: 801-588-0497

Echelon-inc.co