

Echelon Biosciences Inc.

Select-HA LADDERS™



Product Name	Catalog Number	Molecular Mass	Size
Select-HA™ MegaLadder	HYA-MGLAD-20	2 MDa – 8 MDa*	25 µg
Select-HA™ HiLadder	HYA-HILAD-20	500 kDa – 1500 kDa*	25 µg
Select-HA™ LoLadder	HYA-LOLAD-20	30 kDa – 500 kDa*	25 µg

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 synthase¹. Select-HA™ HiLadder and LoLadder contain 5 Select-HA™ molecular mass markers while the Select-HA™ MegaLadder is a mixture of four biotin Select-HA complexes. Select-HA™ is a trademark of Hyalose LLC.

Properties:

Size – 20 lanes. See corresponding amount in table above.

Form – HiLadder and LoLadder: lyophilized (vial may appear empty), MegaLadder: solution of hyaluronan polymer sodium salts

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

Reconstitution – Centrifuge the tube for a few seconds to collect the Select-HA™ solids in the bottom of the tube. Carefully open and add 100 µL of sterile water directly to the bottom of the tube. Allow two hours at 4 °C for sample rehydration and then mix well before use. No reconstitution is required for the MegaLadder.

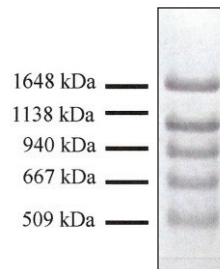
Other – Low voltage (30-40 V) agarose gel runs are recommended for running the Select-HA™ HiLadder and MegaLadder to insure clearly defined bands.

*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. Various biological activities are influenced by the HA size or chain length including proliferation, angiogenesis, inflammation, and binding.

Data: Agarose Gel



Electrophoresis of 5 µL reconstituted Select-HA Ladder™ using a standard gel loading buffer on an agarose gel (0.6 – 1%) results in clearly defined bands when stained with 0.005% Stains-all (in 50% ethanol)². Please see Certificate of Analysis for lot specific ladder M.W.

References:

1. Jing, W.; DeAngelis, P. L. (2004) Synchronized chemoenzymatic synthesis of monodisperse hyaluronan polymers. *J Biol Chem*, 279 (40), 42345-9.
2. Lee, H. G.; Cowman, M. K. (1994) An agarose gel electrophoretic method for analysis of hyaluronan molecular weight distribution. *Anal Biochem*, 219 (2), 278-87.

Related Products:

Product	Catalog Number
Compounds	
BODIPY-HA Texas Red-HA Select-HA™	H-025F, H-250F, H-700F H-025R, H-250R, H-700R HYA-50kEF-1, HYA-500kEF-1, HYA-601kEF-1, HYA-1000kEF-1
Biotinylated Select-HA™	HYA-B50-200, HYA-B250-200, HYA-B500-200, HYA-B1000-200
nanoHA™ HAase Inhibitor	HYA-NAN05-1 B-0601
HA Binding Proteins	
Versican G1 Domain	G-HA01, G-HA02
Assays	
HA Quantification ELISAs Hyaluronidase Activity ELISA	K-1200, K-4800 K-6000

Technical Data Sheet Rev. 3, 10-27-20 - 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.com