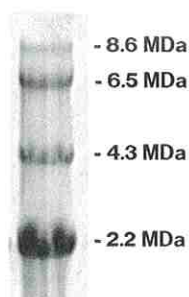


# Certificate of Analysis

<b>Product Name:</b>	Select-HA Mega Ladder				
<b>Product Number:</b>	HYA-MGLAD				
<b>Lot Number:</b>	FVC071824				
<b>Product Description:</b>	0.1 mL solution containing 4 biotin-HA/streptavidin complex with molecular weight range between approximately 2 MDa and 8 MDa				
<b>Quality Test:</b>	<ul style="list-style-type: none"> <li>• Electrophoresis</li> <li>• Multi-angle Laser Light Scattering-Size Exclusion Chromatography (MALLS-SEC)</li> </ul>				
<b>Quality Parameters:</b>	<u>Parameter</u>		<u>Actual</u>	<u>Required</u>	<u>Pass</u>
	A. Markers Clarity on Gel		Clear	Clear	✓
	B. Polydispersity				
	Mn	Mw	Mw/Mn	Required	Pass
	2141 kDa	2151 kDa	1.005	≤1.05	✓
	<p><b>Figure 1. Electrophoresis Analysis</b>            5 µL of Select-HA Mega Ladder (Lot# FVC071824) was run on 0.5% (w/v) TAE agarose gel and stained with 0.005% Stains-All in 50% ethanol</p>				

**This product is for research use only. Not intended or approved for diagnostic or therapeutic use.**

Passes All Functional Specifications:  Angie M. Branch, Biology Manager

Date: 7/26/2024

**For research use only. Not intended for diagnostic or therapeutic use.**