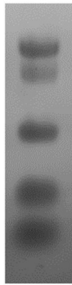
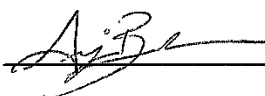


# Echelon Biosciences Inc.

## Certificate of Analysis

<b>Product Name:</b>	Select-HA LoLadder																																			
<b>Product Number:</b>	HYA-LOLAD																																			
<b>Lot Number:</b>	FVC010422																																			
<b>Product Description:</b>	Lyophilized white solids containing 5 Select-HA markers with molecular weight range from 30 kDa to 500 kDa																																			
<b>Quality Assay:</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>	<table border="1"><thead><tr><th>Parameter</th><th>Required</th><th>Actual</th><th>Pass</th></tr></thead><tbody><tr><td>A. Markers Clarity on Gel</td><td>Clear</td><td>Clear</td><td>✓</td></tr><tr><td>B. Polydispersity</td><td></td><td></td><td></td></tr><tr><td>Marker #1</td><td>≤1.05</td><td>1.025</td><td>✓</td></tr><tr><td>Marker #2</td><td>≤1.05</td><td>1.010</td><td>✓</td></tr><tr><td>Marker #3</td><td>≤1.05</td><td>1.011</td><td>✓</td></tr><tr><td>Marker #4</td><td>≤1.05</td><td>1.005</td><td>✓</td></tr><tr><td>Marker #5</td><td>≤1.05</td><td>1.014</td><td>✓</td></tr></tbody></table>	Parameter	Required	Actual	Pass	A. Markers Clarity on Gel	Clear	Clear	✓	B. Polydispersity				Marker #1	≤1.05	1.025	✓	Marker #2	≤1.05	1.010	✓	Marker #3	≤1.05	1.011	✓	Marker #4	≤1.05	1.005	✓	Marker #5	≤1.05	1.014	✓	<p><b>Figure 1. Electrophoresis Analysis</b> 5 µL of reconstituted Select-HA LoLadder (Lot# FVC010422; 1 vial reconstituted in 100 µL water) was run on 1% (w/v) agarose gel and stained with 0.005% Stains-all in 50% ethanol.</p>  <p>Marker #1 – 480 kDa Marker #2 – 400 kDa Marker #3 – 198 kDa Marker #4 – 74 kDa Marker #5 – 32 kDa</p>		
Parameter	Required	Actual	Pass																																	
A. Markers Clarity on Gel	Clear	Clear	✓																																	
B. Polydispersity																																				
Marker #1	≤1.05	1.025	✓																																	
Marker #2	≤1.05	1.010	✓																																	
Marker #3	≤1.05	1.011	✓																																	
Marker #4	≤1.05	1.005	✓																																	
Marker #5	≤1.05	1.014	✓																																	

Passes All Functional Specifications:



Date: 01/19/2022

Angie M. Branch, Biology Manager

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

Ph: 866-588-0455  
Fax: 801-588-0497

Echelon-inc.com