


Echelon Biosciences Inc.

Certificate of Analysis

Product Name:	Select-HA HiLadder																																																					
Product Number:	HYA-HiLAD																																																					
Lot Number:	FVC011023A																																																					
Product Description:	Lyophilized white solids containing 5 Select-HA markers with molecular weight range between approximately 500 kDa and 1500 kDa																																																					
Quality Assay:	<ul style="list-style-type: none"> • Electrophoresis • Multi-angle Laser Light Scattering-Size Exclusion Chromatography (MALLS-SEC) 																																																					
Quality Parameters:	<table border="1"> <thead> <tr> <th>Parameter</th> <th>Actual</th> <th>Required</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 colspan="4">B. Polydispersity</td> </tr> <tr> <th>Marker</th> <th>Mn</th> <th>Mw</th> <th>Mw/Mn</th> <th>Required</th> <th>Pass</th> </tr> <tr> <td>#1</td> <td>1271 kDa</td> <td>1272 kDa</td> <td>1.001</td> <td>≤1.05</td> <td>✓</td> </tr> <tr> <td>#2</td> <td>932 kDa</td> <td>932 kDa</td> <td>1.000</td> <td>≤1.05</td> <td>✓</td> </tr> <tr> <td>#3</td> <td>762 kDa</td> <td>763 kDa</td> <td>1.001</td> <td>≤1.05</td> <td>✓</td> </tr> <tr> <td>#4</td> <td>547 kDa</td> <td>547 kDa</td> <td>1.000</td> <td>≤1.05</td> <td>✓</td> </tr> <tr> <td>#5</td> <td>468 kDa</td> <td>480 kDa</td> <td>1.026</td> <td>≤1.05</td> <td>✓</td> </tr> </tbody> </table> <div style="display: flex; align-items: flex-start;"> <div style="flex: 1;"> <p>Figure 1. Electrophoresis Analysis 5 µL of reconstituted Select-HA HiLadder (Lot# FVC011023A; 1 vial reconstituted in 100 µL water) was run on 0.75% (w/v) agarose gel and stained with 0.005% Stains-all in 50% ethanol.</p> </div> <div style="flex: 0.5; text-align: center;">  </div> <div style="flex: 1;"> <p>Marker #1 – 1272 kDa Marker #2 – 932 kDa Marker #3 – 762 kDa Marker #4 – 547 kDa Marker #5 – 480 kDa</p> </div> </div>						Parameter	Actual	Required	Pass	A. Markers Clarity on Gel	Clear	Clear	✓	B. Polydispersity				Marker	Mn	Mw	Mw/Mn	Required	Pass	#1	1271 kDa	1272 kDa	1.001	≤1.05	✓	#2	932 kDa	932 kDa	1.000	≤1.05	✓	#3	762 kDa	763 kDa	1.001	≤1.05	✓	#4	547 kDa	547 kDa	1.000	≤1.05	✓	#5	468 kDa	480 kDa	1.026	≤1.05	✓
Parameter	Actual	Required	Pass																																																			
A. Markers Clarity on Gel	Clear	Clear	✓																																																			
B. Polydispersity																																																						
Marker	Mn	Mw	Mw/Mn	Required	Pass																																																	
#1	1271 kDa	1272 kDa	1.001	≤1.05	✓																																																	
#2	932 kDa	932 kDa	1.000	≤1.05	✓																																																	
#3	762 kDa	763 kDa	1.001	≤1.05	✓																																																	
#4	547 kDa	547 kDa	1.000	≤1.05	✓																																																	
#5	468 kDa	480 kDa	1.026	≤1.05	✓																																																	

Passes All Functional Specifications: AB

Date: 02/08/2023

Angie M. Branch, Biology Manager

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