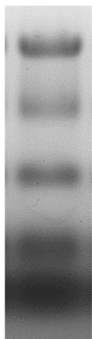
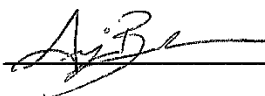


Echelon Biosciences Inc.

Certificate of Analysis

Product Name:	Select-HA LoLadder																																			
Product Number:	HYA-LOLAD																																			
Lot Number:	XCM052824																																			
Product Description:	Lyophilized white solids containing 5 Select-HA markers with molecular weight range from 30 kDa to 550 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>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>1.00</td><td>✓</td></tr><tr><td> Marker #2</td><td>≤1.05</td><td>1.01</td><td>✓</td></tr><tr><td> Marker #3</td><td>≤1.05</td><td>1.01</td><td>✓</td></tr><tr><td> Marker #4</td><td>≤1.05</td><td>1.01</td><td>✓</td></tr><tr><td> Marker #5</td><td>≤1.05</td><td>1.01</td><td>✓</td></tr></tbody></table>	Parameter	Required	Actual	Pass	A. Markers Clarity on Gel	Clear	Clear	✓	B. Polydispersity				Marker #1	≤1.05	1.00	✓	Marker #2	≤1.05	1.01	✓	Marker #3	≤1.05	1.01	✓	Marker #4	≤1.05	1.01	✓	Marker #5	≤1.05	1.01	✓			
Parameter	Required	Actual	Pass																																	
A. Markers Clarity on Gel	Clear	Clear	✓																																	
B. Polydispersity																																				
Marker #1	≤1.05	1.00	✓																																	
Marker #2	≤1.05	1.01	✓																																	
Marker #3	≤1.05	1.01	✓																																	
Marker #4	≤1.05	1.01	✓																																	
Marker #5	≤1.05	1.01	✓																																	
	<p>Figure 1. Electrophoresis Analysis</p> <p>5 µL of reconstituted Select-HA LoLadder (Lot# XCM052824; 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 – 547 kDa</p> <p>Marker #2 – 400 kDa</p> <p>Marker #3 – 198 kDa</p> <p>Marker #4 – 74 kDa</p> <p>Marker #5 – 32 kDa</p>																																	

Passes All Functional Specifications:



Date: 06/04/2024

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