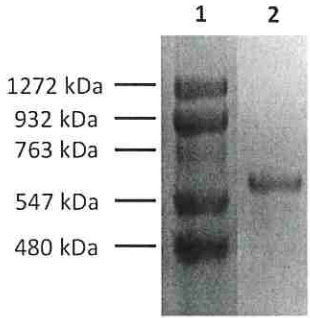


# Echelon Biosciences Inc.

## Certificate of Analysis

<b>Product Name:</b>	600 kDa Select-HA			
<b>Product Number:</b>	HYA-0600			
<b>Lot Number:</b>	FVC031323A			
<b>Product Description:</b>	Lyophilized white solids containing 1 mg Select-HA (determined by Carbazole Assay)			
<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> <li>• Chromogenic Limulus Amebocyte Lysate (LAL) Assay</li> </ul>			
<b>Quality Parameters</b>	<b>Parameter</b>	<b>Required</b>	<b>Actual</b>	<b>Pass</b>
	A. Clarity on Gel	Clear	Clear	✓
	B. Polydispersity	≤1.05	1.000	✓
	C. Molecular Mass	550-650 kDa	641 kDa	✓
	D. Endotoxin Level	≤0.1 EU/mg	0.06 EU/mg	✓
	 <p>1      2</p> <p>1272 kDa — 932 kDa — 763 kDa — 547 kDa — 480 kDa —</p>	<p><b>Electrophoresis Analysis</b> Select-HA was run on 1% (w/v) agarose gel and stained with 0.005% Stains-All in 50% ethanol.</p> <p><b>Lane 1</b> – Select-HA HiLadder (HYA-HILAD, lot# FVC011023A)</p> <p><b>Lane 2</b> – 600 kDa Select-HA (HYA-0600, lot# FVC031323A)</p>		

Passes All Functional Specifications:



Date: 6/22/23

Angie M. Branch, Biology Manager

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