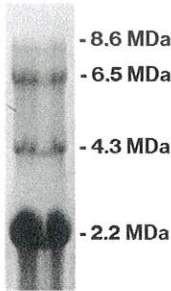


Certificate of Analysis

Product Name:	Select-HA Mega Ladder																											
Product Number:	HYA-MGLAD																											
Lot Number:	FVC081524																											
Product Description:	0.1 mL solution containing 4 biotin-HA/streptavidin complex with molecular weight range between approximately 2 MDa and 8 MDa																											
Quality Test:	<ul style="list-style-type: none"> • Electrophoresis • Multi-angle Laser Light Scattering-Size Exclusion Chromatography (MALLS-SEC) 																											
Quality Parameters:	<table style="width: 100%; border-collapse: collapse;"> <thead> <tr> <th style="text-align: left;">Parameter</th> <th style="text-align: center;">Actual</th> <th style="text-align: center;">Required</th> <th style="text-align: center;">Pass</th> </tr> </thead> <tbody> <tr> <td>A. Markers Clarity on Gel</td> <td style="text-align: center;">Clear</td> <td style="text-align: center;">Clear</td> <td style="text-align: center;">✓</td> </tr> <tr> <td>B. Polydispersity</td> <td colspan="4"></td> </tr> <tr> <td style="padding-left: 20px;">Mn</td> <td style="text-align: center;">Mw</td> <td style="text-align: center;">Mw/Mn</td> <td style="text-align: center;">Required</td> <td style="text-align: center;">Pass</td> </tr> <tr> <td style="padding-left: 20px;">2141 kDa</td> <td style="text-align: center;">2151 kDa</td> <td style="text-align: center;">1.005</td> <td style="text-align: center;">≤1.05</td> <td style="text-align: center;">✓</td> </tr> </tbody> </table>					Parameter	Actual	Required	Pass	A. Markers Clarity on Gel	Clear	Clear	✓	B. Polydispersity					Mn	Mw	Mw/Mn	Required	Pass	2141 kDa	2151 kDa	1.005	≤1.05	✓
	Parameter	Actual	Required	Pass																								
A. Markers Clarity on Gel	Clear	Clear	✓																									
B. Polydispersity																												
Mn	Mw	Mw/Mn	Required	Pass																								
2141 kDa	2151 kDa	1.005	≤1.05	✓																								
Quality Parameters:	<p>Figure 1. Electrophoresis Analysis</p> <p>5 µL of Select-HA Mega Ladder (Lot# FVC081524) 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: 8/23/2024

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