



Certificate of Analysis

Mega-HA™ Ladder

Lot No.: SML200903A

Appearance

Clear liquid (shipped frozen)

Quantity

Vial with 80 µl aqueous solution containing ~5 µg of HA (determined by Carbazole Assay based on free acid form). One vial sufficient for 20 lanes in mini-gel format.

Molecular Masses

Values as determined by Multi-angle Laser Light Scattering-Size Exclusion Chromatography (MALLS-SEC) are typically within $\pm 5\%$. Results here are the average of duplicate analysis.

Mn, number-average molecular mass;

Mw, weight-average molecular mass;

Mw/Mn, polydispersity.

Mn	Mw	Mw/Mn
1949 k	2025 k	1.067

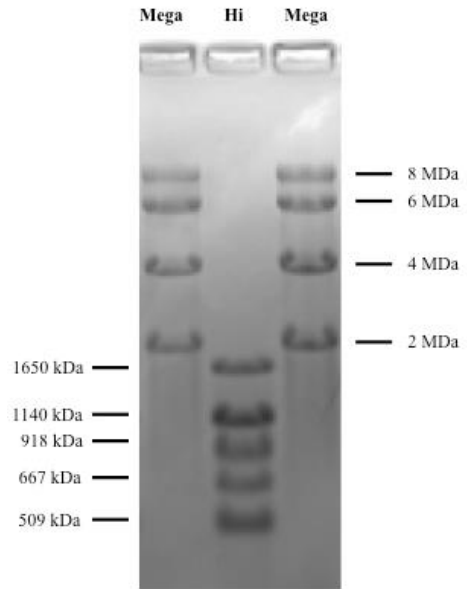


Figure 1. Electrophoretic analysis of Select-HA Mega & Hi-Ladder. Four µl of the thawed Mega-HA™ Ladder solution (Lot# 200903A) or five µl of reconstituted Select-HA™ HiLadder (Lot HL201812; 1 vial in 100 µl total) was run on 0.5% (w/v) agarose gel (mini-gel format) and stained with Stains-all (0.005% Stains-all in 50% ethanol overnight with gentle shaking, destained with water for 2 days). **Mw** molecular masses are indicated.

QC Acceptance Date

Aug 2019

Paul DeAngelis, Ph.D.

Dixy Green, B.S.