

## Hyaluronic Acid (HA)

Echelon Biosciences' HA assay product line offers different assays for HA detection. Refer to the table below to find the assay best fits your needs. See our website for other related products, assay service, related assays.

Support: [echelon@echelon-inc.com](mailto:echelon@echelon-inc.com)

Assay Format	Competitive ELISA	Sandwich ELISA	AlphaScreen™
Catalog Number	K-1200	K-4800	K-5800
No. of wells/kit	96	96	500
Applications	Measures HA concentration in biological samples	Measures HA concentration in biological samples	Measures purified HA concentration from biological sources
Assay Running Time	2.5 – 3 hours	2.5 – 3 hours	1.5 hours
Incubation Temperature	4 °C, 37 °C & Room Temperature	Room Temperature	Room Temperature
HA Standards	Purified HA with mixed HA sizes	Purified HA with mixed HA sizes	Purified HA with mixed HA sizes
Size of HA that is Detected	All Sizes	>130 kDa	All Sizes
HA Detection Range	50-1600 ng/mL	12.5-3200 ng/mL	1.56-1600 ng/mL
Lower Limit of Detection	25 ng/mL	25 ng/mL	15 ng/mL
Sample Volume (duplicate)	100 – 200 µL	25 µL	10 µL
Intra & Inter Assay CVs	<5% Intra CV <10% Inter CV	<5% Intra CV <20% Inter CV	N/A