

Hyaluronic Acid (HA) Assay Comparison

Description:

A noninvasive assessment marker for liver & rheumatoid disease.

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, etc.

Support: 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 (Glycobiology. 2011Feb;21(2):175-83.)	>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	40 ng/mL	5 ng/mL	20 ng/mL
Sample Volume (duplicate)	100– 200µL	25 µL	10µL
Intra & Inter Assay CVs	<20% Intra CV <10% Inter CV (Glycobiology. 2011Feb;21(2):175-83.)	<5% Intra CV <20% Inter CV	N/A

*LoD is defined as the lowest concentration with 2 standard deviations greater than the Limit of Blank (LoB). LoD is calculated by $LoD = LoB + 1.645(SD_{low\ concentration\ sample})$. LoB is calculated by $LoB = mean_{blank} + 1.645(SD_{blank})$.