

# Echelon Biosciences Inc.

## Hyaluronic Acid (HA)

### 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,

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<br><i>(Glycobiology.2011 Feb;21(2):175-83.)</i>                      | >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     | <20% Intra CV<br><10% Inter CV<br><i>(Glycobiology.2011 Feb;21(2):175-83.)</i> | <5% Intra CV<br><20% Inter CV                   | N/A  |

Technical Data Sheet Rev. 4, 01-18-21 - For research use only. Not intended or approved for diagnostic or therapeutic



Echelon Biosciences Inc.

Ph: 866-588-0455

Fax: 801-588-0497

[Echelon-inc.com](http://Echelon-inc.com)