

Simoa® HD-1/HD-X Analyzer

Ultrasensitive Biomarker Assay Menu

Singleplex Advantage Assays - Human

| Analyte | LoD (pg/mL) | LoQ (pg/mL) | Dynamic Range (pg/mL) | Median Endogenous (pg/mL) | Sample Volume† | Sample Type†† | Catalog Number |
|---|-------------|-------------|-----------------------|--|----------------|---------------|----------------|
| Aβ40 | 0.522 | 1.23 | 0-800 | 65.97 | 32.5 µL | C, E | 101672 |
| Aβ42 | 0.044 | 0.137 | 0-400 | 4.7 | 32.5 µL | C, E | 101664 |
| C-Peptide | 0.013 | 0.021 | 0-400 | 1559 | 32.5 µL | E, S | 100199 |
| Eotaxin/CCL-11 | 0.1 | 0.18 | 0-960 | 75.6 | 32.5 µL | E, S | 101212 |
| G-CSF | 0.095 | 0.095 | 0-400 | 7.12 | 32.5 µL | E, S | 101235 |
| GM-CSF | 0.0019 | 0.0103 | 0-120 | 0.0865 | 32.5 µL | E, S | 102329 |
| IFNα | 0.0025 | 0.0047 | 0-150 | 0.0036 | 100 µL | S | 100860 |
| IFNγ | 0.0155 | 0.098 | 0-100 | 0.522 | 65 µL | E, S | 103337 |
| IL-1β | 0.016 | 0.083 | 0-240 | 0.058 | 100 µL | E, S | 101605 |
| IL-2 | 0.011 | 0.041 | 0-120 | 0.086 | 49.5 µL | E, S | 101635 |
| IL-4 | 0.0046 | 0.039 | 0-200 | 0.024 | 65 µL | E, S | 100196 |
| IL-5 | 0.004 | 0.0165 | 0-12 | 0.22 | 91 µL | E, S | 102860 |
| IL-6 | 0.0055 | 0.01 | 0-120 | 1.73 | 32.5 µL | E, S | 101622 |
| IL-7 | 0.009 | 0.103 | 0-600 | 28.0 | 32.5 µL | E, S | 103277 |
| IL-8 | 0.056 | 0.0921 | 0-1,200 | 5.31 | 32.5 µL | E, S | 100198 |
| IL-10 | 0.0038 | 0.021 | 0-120 | 0.94 | 32.5 µL | E, S | 101643 |
| IL-12p70 | 0.0048 | 0.017 | 0-40 | 1.95 | 32.5 µL | E, S | 100988 |
| IL-13 | 0.002 | 0.005 | 0-30 | 0.039 | 65 µL | E, S | 102732 |
| IL-15 | 0.003 | 0.0062 | 0-40 | 3.23 | 32.5 µL | E, S | 100794 |
| IL-17A | 0.0042 | 0.021 | 0-120 | 0.124 | 32.5 µL | E, S | 101599 |
| IP-10 | 0.052 | 0.177 | 0-800 | 105 | 32.5 µL | E, S | 101132 |
| MCP-1 | N/A | 0.153 | 0-800 | 85.3 | 32.5 µL | E, S | 101154 |
| NF-light™ | 0.038 | 0.174 | 0-2,000 | 5.33 | 45.5 µL | C, E, S | 102258 |
| HIV p24 | 0.0027 | 0.01 | N/A | N/A | 154 µL | E, S | 102215 |
| PSA | 0.015 | 0.024 | 0-400 | 1.81 | 32.5 µL | E, S | 101478 |
| Tau | 0.019 | 0.061 | 0-360 | 1.65 | 45.5 µL | C, E, S | 101552 |
| SARS-CoV-2 Spike IgG (HD-X Only) | 15.0 | 8,200 | 0-280,000,000 | 32,000 (pre-pand) 3,360,000 (PCR Pos) | 5 µL | E, S | 103769 |
| SARS-CoV-2 N Protein Antigen (HD-X Only) | 0.099 | 0.313 | 0-800 | 2,338 (PCR Pos) | 32.5 µL | NP Swab | 103806 |
| SNAP-25 | 0.880 | 2.56 | 0-1,000 | 59.6 | 32.5 µL | C | 103575 |
| pTau-181 | 0.724 | 2.640 | 0-85 | 32.0 | 32.5 µL | C | 103377 |
| pTau-181 V2 | 0.028 | 0.085 | 0-412 | 0.856 | 32.5 µL | C, E, S | 103714 |
| pTau-231 | 0.621 | 1.83 | 0-1,200 | 20.8 | 45.5 µL | C | 102292 |
| TDP-43 | 2.48 | 8.23 | 0-8,000 | 130 | 32.5 µL | C, E, S | 103293 |
| TNF-α | 0.016 | 0.034 | 0-400 | 1.94 | 32.5 µL | E, S | 101580 |
| TRAIL | 0.0083 | 0.0177 | 0-400 | 23.1 | 32.5 µL | E, S | 100906 |

† Volumes shown are for at bench dilutions. On-instrument dilutions require higher sample volumes | †† C = CSF, E = EDTA plasma, S = serum

Multiplex Advantage Assays - Human

| Assay | Analytes | LoD (pg/mL) | LoQ (pg/mL) | Dynamic Range (pg/mL) | Median Endogenous (pg/mL) | Sample Volume [†] | Sample Type ^{**} | Catalog Number |
|---------------------------|----------|-------------|-------------|-----------------------|---------------------------|----------------------------|---------------------------|----------------|
| Neurology 2-Plex B | GFAP | 0.475 | 4.15 | 0-40,000 | 36.9 | 32.5 µL 5 µL (CSF) | C, E, S | 103520 |
| | Nf-L | 0.065 | 0.20 | 0-2000 | 3.63 | | | |
| Neurology 3-Plex A | Aβ40 | 0.196 | 0.675 | 0-800 | 209 | 45.5 µL 5 µL (CSF) | C, E | 101995 |
| | Aβ42 | 0.045 | 0.142 | 0-400 | 11.1 | | | |
| | Tau | 0.019 | 0.063 | 0-400 | 1.43 | | | |
| Neurology 4-Plex A | GFAP | 0.221 | 0.467 | 0-4,000 | 89.7 | 45.5 µL 5 µL (CSF) | C, E, S | 102153 |
| | Nf-L | 0.104 | 0.241 | 0-2000 | 10.6 | | | |
| | Tau | 0.024 | 0.053 | 0-400 | 2.21 | | | |
| | UCH-L1 | 1.74 | 5.45 | 0-40,000 | 12.21 | | | |
| Neurology 4-Plex B | GFAP | 1.18 | 9.38 | 0-40,000 | 50.0 | 32.5 µL 5 µL (CSF) | C, E, S | 103345 |
| | Nf-L | 0.096 | 0.500 | 0-2,000 | 5.91 | | | |
| | Tau | 0.037 | 0.125 | 0-400 | 0.95 | | | |
| | UCH-L1 | 2.43 | 9.38 | 0-40,000 | 6.81 | | | |
| Neurology 4-Plex E | Aβ40 | 0.384 | 1.02 | 0-280 | 89.4 | 32.5 µL 5 µL (CSF) | C, E | 103670 |
| | Aβ42 | 0.136 | 0.378 | 0-100 | 5.69 | | | |
| | GFAP | 0.441 | 2.89 | 0-20,000 | 52.2 | | | |
| | Nf-L | 0.090 | 0.400 | 0-2,000 | 10.0 | | | |

[†] Volumes shown are for at bench dilutions. On-instrument dilutions require higher sample volumes | ^{**} C = CSF, E = EDTA plasma, S = serum

Simoa Discovery assay kits are provided as complete kits packaged with reagents for processing up to 192 samples. Includes calibrator concentrate, calibrator diluent, bead stock, bead diluent, detector stock, detector diluent, SBG concentrate, SBG diluent, RGP, and sample diluent. Representative performance is determined in biological matrices and includes:

- Dynamic range
- Sensitivity
- Calibration curve fitting
- Intra-run, inter-plate, and inter-lot precision and accuracy
- Robustness and stability of calibrators, antibodies, and controls
- Analysis of normal samples for the specified species to determine the normal range of biomarker concentration detected with the assay
- Specificity, spike recovery, and dilution linearity

Discovery Assays – Human

| Analyte | LoD (pg/mL) | LoQ (pg/mL) | Dynamic Range (pg/mL) | Median Endogenous (pg/mL) | Sample Volume [†] | Sample Type ^{††} | Catalog Number |
|----------------------|-------------|-------------|-----------------------|---------------------------|----------------------------|---------------------------|----------------|
| α-Synuclein | 0.955 | 4.12 | 0-10,000 | 4,145 | 18.2 µL | C, E, S | 102233 |
| BDNF | 0.011 | 0.034 | 0-64,000 | 11,306 | 5 µL | C, E, S | 102039 |
| CA-125 | 0.003 U/mL | 0.010 U/mL | 0-200 U/mL | 1.62 U/mL | 13 µL | E, S | 102136 |
| Cathepsin S | 0.7 | 1.95 | 0-200 | 6,566 | 5 µL | E, S | 102064 |
| CEA | 0.486 | 2.33 | 0-85 ng/mL | 1511.5 | 5 µL | E, S | 102556 |
| c-MET | 0.036 | 0.244 | 0-800 | 58,012 | 5 µL | E, L, S | 102073 |
| CRP | 0.048 | 0.686 | 0-48 | 19 | 5 µL | E, S | 102583 |
| CXCL13 | 0.048 | 0.07 | 0-800 | 22.63 | 32.5 µL | E, S | 102635 |
| GFAP | 0.211 | 0.686 | 0-4,000 | 88 | 45.5 µL | C, E, S | 102336 |
| HE4 / WFDC2 | 0.135 | 0.977 | 0-4,000 | 104 | 32.5 µL | E, S | 103059 |
| IL-1α | 0.004 | 0.01 | 0-60 | 0.0293 | 65 µL | E, S | 101968 |
| IL-3 | 0.226 | 0.686 | 0-2000 | 0.279 | 45.5 µL | E, S | 102462 |
| IL-17C | 0.065 | 0.206 | 0-1,200 | 1.66 | 32.5 µL | E, S | 102570 |
| IL-18 | 0.004 | 0.012 | 0-2250 | 200.1 | 5 µL | C, E, S | 102700 |
| IL-22 (Total) | 0.0054 | 0.0103 | 0-120 | 7.16 | 32.5 µL | E, S | 103071 |
| IL-28A | 0.022 | 0.069 | 0-200 | 0.303 | 65 µL | E, S | 101419 |
| IL-36β | 0.01 | 0.206 | 0-600 | 0.426 | 32.5 µL | E, S | 101808 |
| Leptin | 2.46 | 4.94 | 0-60,000 | 6,083 | 5 µL | E, S | 101855 |
| LIF | 0.015 | 0.086 | 0-520 | 0.412 | 32.5 µL | E, S | 102394 |
| MCP-3 | 0.124 | 0.309 | 0-450 | 0.445 | 75 µL | E, S | 102382 |
| MMP-9 | 0.581 | 4.88 | 0-5,000 ng/ml | 555 | 5 µL | C, E, S | 102491 |
| pNF-heavy | 0.663 | 2.88 | 0-8,400 | 30.82 | 32.5 µL | C, E, S | 102669 |
| NSE | 1.296 | 9.88 | 0-120 | 7,845 | 5 µL | C, E, S | 102475 |
| PD-1 | 0.247 | 0.879 | 0-7200 | 73 | 32.5 µL | E, S | 102929 |
| PD-L1 | 0.055 | 0.617 | 0-4,300 | 33.79 | 6.5 µL | E, S | 102648 |
| PIGF | 0.064 | 0.3 | 0-960 | 3.82 | 37.5 µL | E, S | 102318 |
| TGFα | 0.031 | 0.207 | 0-900 | 3.34 | 65 µL | E, S | 101863 |
| TGFβ | 0.137 | 0.514 | 0-24,000 | 34,836 | 8.5 µL | E, S | 101984 |
| UCH-L1 | 1.05 | 3.43 | 0-20,000 | 9.51 | 45.5 µL | C, E, S | 102343 |
| VEGF | 0.041 | 0.137 | 0-800 | 3.49 | 32.5 µL | E, S | 102794 |

[†] Volumes shown are for at bench dilutions. On-instrument dilutions require higher sample volumes | ^{††} C = CSF, E = EDTA plasma, S = serum

Simoa Discovery assay kits are provided as complete kits packaged with reagents for processing up to 192 samples. Includes calibrator concentrate, calibrator diluent, bead stock, bead diluent, detector stock, detector diluent, SBG concentrate, SBG diluent, RGP, and sample diluent. Representative performance is determined in biological matrices and includes:

- Dynamic range
- Sensitivity
- Calibration curve fitting
- Intra-run, inter-plate, and inter-lot precision and accuracy
- Robustness and stability of calibrators, antibodies, and controls
- Analysis of normal samples for the specified species to determine the normal range of biomarker concentration detected with the assay
- Specificity, spike recovery, and dilution linearity

Discovery Assays – Mouse

| Analyte | LoD (pg/mL) | LoQ (pg/mL) | Dynamic Range (pg/mL) | Median Endogenous (pg/mL) | Sample Volume [†] | Sample Type ^{††} | Catalog Number |
|--------------------------------|-------------|-------------|-----------------------|---------------------------|----------------------------|---------------------------|----------------|
| IL-1α | 0.0085 | 0.060 | 0-700 | 1.063 | 13 μ L | E, S | 102504 |
| IL-1β | 0.004 | 0.021 | 0-120 | 0.201 | 32.5 μ L | E, S | 102517 |
| IL-6 | 0.035 | 0.120 | 0-700 | 16.0 | 32.5 μ L | E, S | 102109 |
| IL-17A | 0.088 | 0.206 | 0-2,400 | 33.76 | 17 μ L | E, S | 101824 |
| IL-17F | 0.102 | 0.412 | 0-2,400 | 11.01 | 32.5 μ L | E, S | 101839 |
| IL-22 | 0.095 | 0.206 | 0-1,200 | 273.3 | 32.5 μ L | E, S | 101919 |
| IL-23 | 0.033 | 0.137 | 0-800 | 4.89 | 32.5 μ L | E, S | 101847 |
| TNF-α | 0.132 | 1.230 | 0-7,200 | 22.6 | 32.5 μ L | E, S | 102100 |
| Tau | 0.615 | 0.823 | 0-2,400 | 26.7 | 32.5 μ L | C, E, S | 102209 |

[†]Volumes shown are for at bench dilutions. On-instrument dilutions require higher sample volumes | ^{††}C = CSF, E = EDTA plasma, S = serum

Simoa Discovery assay kits are provided as complete kits packaged with reagents for processing up to 192 samples. Includes calibrator concentrate, calibrator diluent, bead stock, bead diluent, detector stock, detector diluent, SBG concentrate, SBG diluent, RGP, and sample diluent. Representative performance is determined in biological matrices and includes:

- Dynamic range
- Sensitivity
- Calibration curve fitting
- Intra-run, inter-plate, and inter-lot precision and accuracy
- Robustness and stability of calibrators, antibodies, and controls
- Analysis of normal samples for the specified species to determine the normal range of biomarker concentration detected with the assay
- Specificity, spike recovery, and dilution linearity