

Product Number:	102215
Lot Number:	503682
Expiration:	15-Nov-2023
Platform(s):	HD-1/HD-x & SR-X



Data below represents results generated on the Simoa® HD-X Analyzer.

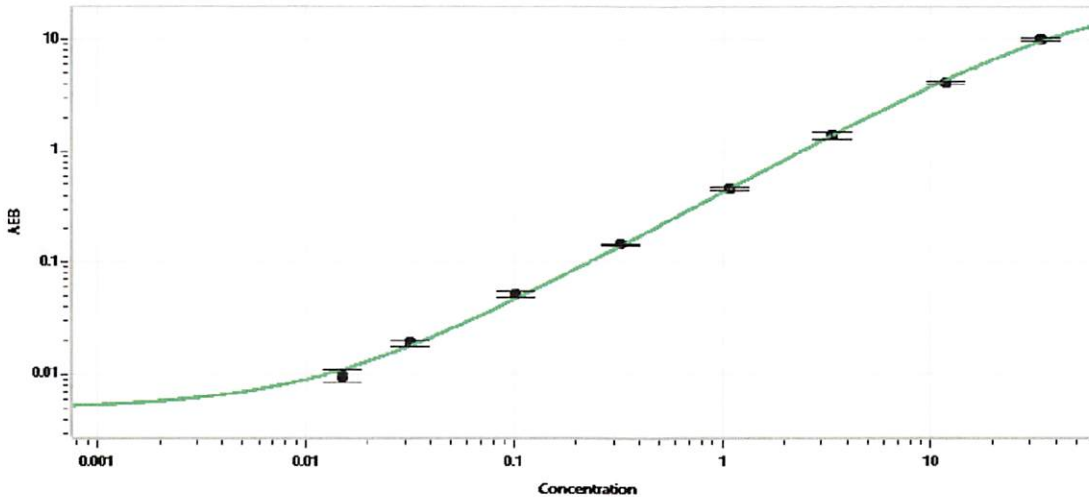
Component	Part Number	Lot Number
Bead Reagent	102211	231912
Detector Reagent	102212	232507
SBG Reagent	102213	232603
Assay Diluent	102214	234809
Calibrators	101310	224201
Control 1	100702	224202
Control 2	100703	224202
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	224202	0.110	0.085-0.127
Control 2	224202	11.6	9.25-13.9
Panel 1	019601	0.204	0.127-0.243
Panel 2	019601	2.04	1.56-2.34
Panel 3	019601	11.7	9.44-14.2

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Review/Approval: Sonia Khillar Name QC Manager Title Sonia Khillar / 21 DEC 2022 Signature/Date