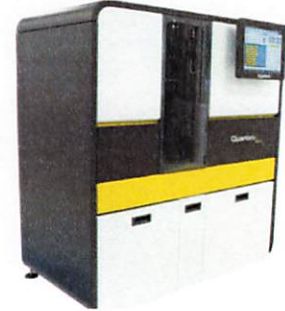


Product Number:	101643
Lot Number:	503891
Expiration:	17-Feb-2024
Platform(s):	HD-1/HD-X and SR-X



Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	101645	315004	315004A
Detector Reagent	101644	315205	315205A
SBG Reagent (HD-1)	101646	315206	315206A
Sample Diluent	102200	308221	N/A
Calibrator Diluent	101648	315304	315304A
Calibrator Concentrate	100670	308004	N/A
SBG Diluent (SR-X) †	103158	N/A ¹	N/A ¹
RGP Reagent	103159	N/A ¹	N/A ¹

¹ These items are not Kit Lot Specific

† Specific for use on the SR-X platform - refer to the Simoa™ Human IL-10 Advantage SR-X Kit Instructions (KI-0036).

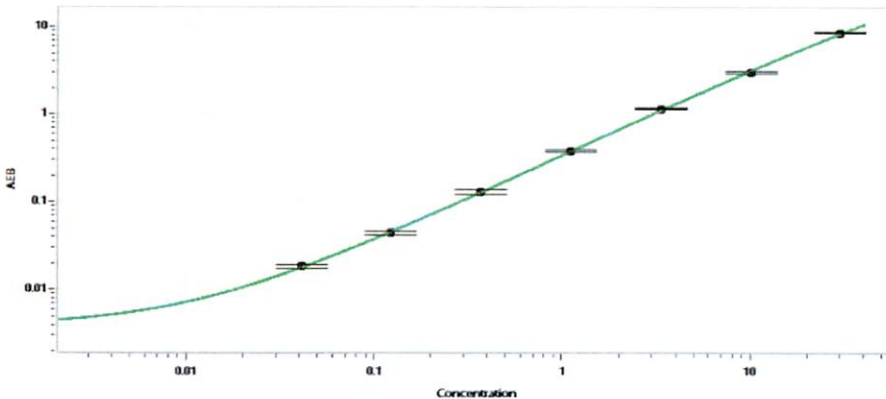
Data below represents results generated on the Simoa® HD-X.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	308004	1.70	1.20 - 2.00
Control 2	308004	24.8	18.0 - 30.0
Panel 1	215202	0.290	0.213 - 0.561
Panel 2	215202	0.536	0.425 - 0.738
Panel 3	306507	53.4	47.1 - 71.5

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

Calibrator Concentrate Value Assignment (pg/mL)	603
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Sonia Khillar
Name

QC
Title

Sonia Ralla 16 JUN 2023
Signature/Date