

Product Number:	104691
Lot Number:	504342
Expiration:	22-Jul-2025
Platform(s):	HD-X



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

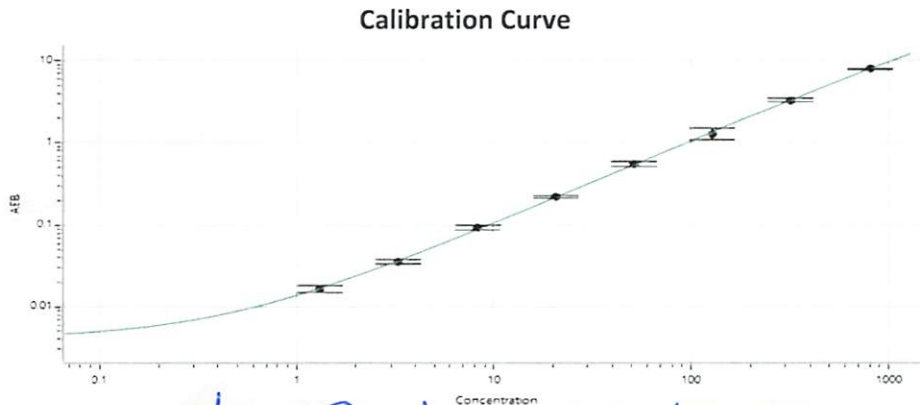
Component	Part Number	Lot Number
Bead Reagent	104752	421812
Detector Reagent	104675	421902
SBG Reagent	104677	421903
Plasma Sample Diluent	104719	421814
CSF Sample Diluent	104720	421815
Calibrator Diluent	104678	402220
Control Diluent	104718	402221
Calibrator Concentrate	104687	402909
Control 1 Concentrate	104689	402910
Control 2 Concentrate	104690	402911
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	402910	137	86.2 - 160
Control 2 Concentrate	402911	1480	1035 - 1922

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

Calibrator Concentrate Reconstitution Volume (uL)	911
Control 1 Concentrate Reconstitution Volume (uL)	911
Control 2 Concentrate Reconstitution Volume (uL)	911



Calibrator Levels (pg/mL)	
A	0.000
B	1.31
C	3.28
D	8.19
E	20.5
F	51.2
G	128
H	320
I	800

Review/Approval

Dhruvi Patel
 Name Title

20 Sep 2024 / DPP
 Signature/Date