

Product Number:	104713
Lot Number:	504371
Paired Reagent Kit Lot Number:	504091
Expiration:	30-Jan-2029
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	104708	403015
Control 1 Concentrate	104710	403016
Control 2 Concentrate	104711	403017

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Review/Approval

Dhruvi Patel Doc Control I QA
 Name Title

DPP / 11 Sep 2024
 Signature/Date