

Product Number:	104628
Lot Number:	504222
Paired Reagent Kit Lot Number:	503964
Expiration:	21-Jun-2028
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	104361	317309
Control 1 Concentrate	104362	317310
Control 2 Concentrate	104363	317311

Calibrator Concentrate Reconstitution Volume (uL)	1244
Control 1 Concentrate Reconstitution Volume (uL)	1244
Control 2 Concentrate Reconstitution Volume (uL)	1244

Review/Approval

ROBERT SOMOGYI QC MANAGER [Signature] 18 APR 2024

---

Name Title Signature/Date