

Product Number:	104692
Lot Number:	504291
Paired Reagent Kit Lot Number:	504093
Expiration:	29-Jan-2029
Platform(s):	HD-X

Component	Part Number	Lot Number
Calibrator Concentrate	104687	402909
Control 1 Concentrate	104689	402910
Control 2 Concentrate	104690	402911



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

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

Review/Approval

Robert Somogyi

QC Manager

RS *24 JUN 2024*

Name

Title

Signature/Date