

Product Number:	104671
Lot Number:	504372
Paired Reagent Kit Lot Number:	504092
Expiration:	29-Jan-2029
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	104666	402913
Control 1 Concentrate	104668	402914
Control 2 Concentrate	104669	402915

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

Review/Approval

Dipak Patil
Name
Doc Control II QA
Title
DP / 10 Sep 2024
Signature/Date