

Product Number:	104571
Lot Number:	504333
Paired Reagent Kit Lot Number:	504307
Expiration:	21-Feb-2029
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	104566	406408
Control 1 Concentrate	104568	406409
Control 2 Concentrate	104569	406410

Calibrator Concentrate Reconstitution Volume (uL)	784
Control 1 Concentrate Reconstitution Volume (uL)	784
Control 2 Concentrate Reconstitution Volume (uL)	784

Review/Approval

Robert Somogyi

QC Manager

*[Handwritten Signature]* 14AUG2024

Name

Title

Signature/Date