

Product Number:	104980
Lot Number:	536659
Paired Reagent Kit Lot Number:	504385
Expiration:	23-Apr-2029
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Calibrator Concentrate	104975	411414
Control 1 Concentrate	104977	411416
Control 2 Concentrate	104978	411417



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

Calibrator Concentrate Reconstitution Volume (uL)	1000
Control 1 Concentrate Reconstitution Volume (uL)	1000
Control 2 Concentrate Reconstitution Volume (uL)	1000

Review/Approval

William Chan

document control specialist I

William Chan 04DEC2025

Name

Title

Signature/Date