

Product Number:	101736
Lot Number:	503516
Expiration:	14-Feb-2023



Data below represents results generated on the Simoa® HD-1 / HD-X Analyzer.

Component	Part Number	Lot Number
RGP Reagent	103159	20450
Means of Analysis		PASS/FAIL
HD-1/HD-X		Pass

Review/Approval

Matt Smart
Name

QC
Title

MA 09 Jun 2022
Signature/Date