

Product Number:	101736
Lot Number:	503471
Expiration:	16-Feb-2023



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

Component	Part Number	Lot Number
RGP Reagent	103159	20470

Means of Analysis	PASS/FAIL
HD-1/HD-X	Pass

Review/Approval

Richa Shrestha

QC Supervisor

[Handwritten Signature] 29 Apr 2023

Name

Title

Signature/Date