

Product Number:	101736
Lot Number:	503695
Expiration:	26-Jul-2023



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

Component	Part Number	Lot Number
RGP Reagent	103159	22070

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

Review/Approval

Matt Smart

QC

MRA 19 Sep 2021

Name

Title

Signature/Date