

Product Number:	101736
Lot Number:	503799
Expiration:	19-Dec-2023



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

Component	Part Number	Lot Number
RGP Reagent	103159	23530

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

Review/Approval

Matt Koubouris

QC Supervisor

[Signature] 06Feb2023

Name

Title

Signature/Date