

Product Number:	101736
Lot Number:	503663
Expiration:	23-May-2023



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

Component	Part Number	Lot Number
RGP Reagent	103159	21430

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

Review/Approval

Matt Smart

Name

QC

Title

MAA

03Aug2022

Signature/Date