

Product Number:	101736
Lot Number:	503543
Expiration:	20-Apr-2023



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

Component	Part Number	Lot Number
RGP Reagent	103159	21100

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

Review/Approval

Sonia Khillan
Name

QC
Title

Sonia Khillan / 13JUL2022
Signature/Date