

Product Number:	101736
Lot Number:	503997
Expiration:	5-Aug-2024



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

Component	Part Number	Lot Number
RGP Reagent	103159	31560

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

Review/Approval

Sonia Khillan

QC Manager

Sonia Khillan 05 OCT 2023

Name

Title

Signature/Date