

Product Number:	101736
Lot Number:	503899
Expiration:	20-May-2024



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

Component	Part Number	Lot Number
RGP Reagent	103159	30790

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

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan 05JUN2023
Signature/Date