

Product Number:	101736
Lot Number:	503946
Expiration:	30-Jul-2024



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

Component	Part Number	Lot Number
RGP Reagent	103159	31500

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

Review/Approval

Pratish SNETI
Name

QC SUPERVISOR
Title

Pratish SNETI 09 AUG 2023
Signature/Date