

Product Number:	101736
Lot Number:	504088
Expiration Date:	30-Dec-24



Component	Part Number	Lot Number
RGP Reagent	103159	33030

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

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

Review/Approval

ROBERT SOMOGYI

QC MANAGER

[Signature] 14 DEC 2023

Name

Title

Signature/Date