

Product Number:	101736
Lot Number:	504149
Expiration Date:	19-Jul-2025



Component	Part Number	Lot Number
RGP Reagent	103159	40190

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

08FEB2024

Name

Title

Signature/Date