

Product Number:	103718
Lot Number:	503897
Expiration:	10-Aug-2023



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

Component	Item Number	Lot Number	Control Number
SBG Concentrate	103650	313007	313007A

SBG Concentrate Concentration (nM)	160
------------------------------------	-----

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Sonia Killian
Name

QC Manager
Title

Sonia Killian 31MA2023
Signature/Date