

Product Number:	101735
Lot Number:	504028
Expiration:	31-Jul-2024



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

Component	Item Number	Lot Number	Control Number
SBG Dilution Buffer	100376	327510	N/A
SBG Concentrate	103397	315713	315713A

Laboratory Analysis

Means of Analysis	Characteristic	Result
HD-1 Analyzer	Functional Test	PASS
SBG Concentrate Concentration (nM)		50

Review/Approval

Sonia Khillan
Name

Senior QC Analyst
Title

Sonia Khillan 31 OCT 2023
Signature/Date