

<b>Product Number:</b>	101735
<b>Lot Number:</b>	503900
<b>Expiration:</b>	28-Mar-2024

Component	Item Number	Lot Number	Control Number
SBG Dilution Buffer	100376	311004	311004A
SBG Concentrate	103397	308711	n/a

**Laboratory Analysis**

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



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

Review/Approval

*Sonia Khillan*  
Name

*QC Manager*  
Title

*Sonia Khillan - 12 JUN 2023*  
Signature/Date