

<b>Product Number:</b>	101735
<b>Lot Number:</b>	503726
<b>Expiration:</b>	26-Jul-2023



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

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	225011
SBG Concentrate	103397	222816

**Laboratory Analysis**

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

**Review/Approval**

*Sonia Khillan*  
Name

*QC Manager*  
Title

*Sonia Khillan / 04 NOV 2022*  
Signature/Date