

<b>Product Number:</b>	103650
<b>Lot Number:</b>	220720
<b>Expiration:</b>	26-Oct-2022
<b>Storage:</b>	2-8°C



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

**Laboratory Analysis**

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	135
------------------------------------	-----

Review/Approval

*Sonia Khillan*

Name

QC

Title

*Sonia Khillan / 10 AUG 2022*

Signature/Date