

<b>Product Number:</b>	103650
<b>Lot Number:</b>	227217
<b>Expiration:</b>	29-Nov-2022
<b>Storage:</b>	2-8 °C



*Data below represents results generated on the Simoa<sup>®</sup> HD-1 Analyzer.*

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>
HD-1 Analyzer	PASS
SBG Concentrate Concentration (nM)	160

Review/Approval

*Sonia Khillan*

Name

*Ac Manager*

Title

*Sonia/20/22 / 13OCT2022*

Signature/Date