

<b>Product Number:</b>	103650
<b>Lot Number:</b>	210115
<b>Expiration:</b>	11-Jul-2022
<b>Storage:</b>	2-8 °C



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

**Laboratory Analysis**

<b>Characteristic</b>	<b>Result</b>
HD-1 Analyzer	PASS

<b>SBG Concentrate Concentration (nM)</b>	134.6
---	-------

Review/Approval

*Matt Smart*  
Name

QC  
Title

*MZA 19 May 2022*  
Signature/Date