

<b>Product Number:</b>	103397
<b>Lot Number:</b>	107422
<b>Expiration:</b>	15-Mar-2022
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

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

<b>SBG Concentrate Concentration (nM)</b>	50
---	----

Review/Approval Shivani Goel QC Manager Shivani Goel 03May2021  
Name Title Signature/Date