

<b>Product Number:</b>	103397
<b>Lot Number:</b>	921713
<b>Expiration:</b>	5-Aug-2020
<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**

Patrizia Stadler

Assoc. Director, QC

*p. Stadler*

21 Feb 2020

**Name**

**Title**

**Signature/Date**