

Product Number:	103397
Lot Number:	108324
Expiration:	24-Mar-2022
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	50
------------------------------------	----

Review/Approval

Shivani Goel  
Name

QC  
Title

Shivani Goel 07 Jun 2021  
Signature/Date