

Product Number:	103397
Lot Number:	101913
Expiration:	19-Jan-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Shivani Goel
Junpil Park

QC Manager
~~QC Supervisor~~

Shivani Goel 26 Feb 2021

Name

SG

Title

SG 26 Feb 2021

Signature/Date