

Product Number:	103397
Lot Number:	101116
Expiration:	11-Jan-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Junpil Park
Name

Supervisor, QC
Title

 03 Feb 2021
Signature/Date