

Product Number:	103650
Lot Number:	129811
Expiration:	25-Jan-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	136
---	-----

Review/Approval

Jungil Park
Name

QC Supervisor
Title

 *17 Nov 2021*
Signature/Date