

Product Number:	103397
Lot Number:	926108
Expiration:	18-Sep-2020
Storage:	2-8°C

Laboratory Analysis

Means of Analysis	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	142.6
---	-------

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler

13 NOV 2019

Signature/Date