

Product Number:	101735
Lot Number:	502444
Expiration:	2-May-2021

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	012309
SBG Concentrate	103397	013908

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

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 25Jun2020
Signature/Date