

Product Number:	101735
Lot Number:	501861
Expiration:	15-Apr-2020

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	916108
SBG Concentrate	103397	910511

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

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

18 JUN 2019

Name

Title

Signature/Date