

Product Number:	101735
Lot Number:	501925
Expiration:	22-Apr-2020

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	919202
SBG Concentrate	103397	911210

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

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler

19 Jul 2019

Signature/Date