

Product Number:	101735
Lot Number:	502552
Expiration:	15-Jul-2021

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	026615
SBG Concentrate	103397	019712

Laboratory Analysis

Means of Analysis	Characteristic	Result
HD-1 Analyzer	Functional Test	PASS
SBG Concentrate Concentration (nM)		50

Review/Approval

Patrizia Stadler	Associate Director, QC	<i>P. Stadler</i>	02 OCT 2020
Name	Title	Signature/Date	