

Product Number:	103297
Lot Number:	100810
Expiration:	8-Jan-2022
Storage:	2-8°C

SBG Stock (nM)	10nM
----------------	------

Review/Approval

Junpil Park

QC, Supervisor



14 Jan 2024

Name

Title

Signature/Date