

Product Number:	103591
Lot Number:	504210
Expiration:	10-Nov-2025

Component	Part Number	Lot Number
Bead Reagent	103581	415201
Detector Reagent	103583	415202
SBG Reagent	103585	415203
Control 1	103589	415603
Control 2	103590	415603
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

Bead Count (beads/mL)	1.99 x 10 ⁷
Bead % Monomeric	93%
Isingle	121

Review/Approval

Cody Carlson
Name

QC Manager
Title


Signature/Date