

<b>Product Number:</b>	104827
<b>Lot Number:</b>	504297
<b>Expiration:</b>	11-Oct-2025

Component	Part Number	Lot Number
Bead Diluent Reagent	104834	410110
Detector Diluent Reagent	104850	412719
SBG Diluent Reagent	104829	410212
Control 1 Diluent Reagent	104841	413409
Control 2 Diluent Reagent	104846	413410
Beads	104833	413809
Detector	104849	413810
SBG	104831	413811
Control 1	104839	413812
Control 2	104844	413813
RGP 2.0 Reagent	104321	N/A <sup>1</sup>
Activation Buffer Reagent	104324	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

Review/Approval

Cody Carlson

QC Manager

Name

Title

 28 Jun 2024  
Signature/Date