

Product Number:	104827
Lot Number:	504627
Expiration:	28-May-2027

Component	Part Number	Lot Number
Bead Diluent Reagent	104834	514811
Detector Diluent Reagent	104850	514813
SBG Diluent Reagent	104829	514816
Control 1 Diluent Reagent	104841	514814
Control 2 Diluent Reagent	104846	514815
Beads	104833	512905
Detector	104849	408720
SBG	104831	407817
Control 1	104839	408722
Control 2	104844	408723
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

Review/Approval

Dipak Patel Sr. Doc. Control-QA at 26 Jun 2025
 Name Title Signature/Date