

Product Number:	104827
Lot Number:	504639
Expiration:	16-Jun-2027

Component	Part Number	Lot Number
Bead Diluent Reagent	104834	516726
Detector Diluent Reagent	104850	516725
SBG Diluent Reagent	104829	516724
Control 1 Diluent Reagent	104841	516728
Control 2 Diluent Reagent	104846	516727
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

Dilip Patel

Name

Sr. Doc. Control - QA

Title

CT 05 AUG 2025

Signature/Date