

**VALIDATION AND AUTOMATION OF PROCESSING THE BODE BUCCAL DNA COLLECTOR USING  
POWERPLEX® 16 HS**

Lana Ramos

Bode Technology Group, 10430 Furnace Rd, Suite 107, Woodridge, VA 22193

PowerPlex® 16 HS designed for direct-to-Amplification reagent designed for more efficient processing of reference samples. The PowerPlex® 16 HS system was launched to support FTA collection devices. At Bode, we have developed an automated process for the utilizing PowerPlex® 16 HS with the Bode Buccal DNA Collectors. During this presentation, we will present on the reasons for a modified protocol to process the collectors, the development and optimization of the protocol, automating the protocol and the results of the validation studies.