

IMPLEMENTING THE DNA IQ™ CASEWORK PRO KIT FOR MAXWELL® 16

Ashley Clark

Johnson County Crime Lab, 6000 Lamar Ave, Mission, KS 66202

The DNA IQ™ Casework Sample Kit, by Promega, was recently updated and changes were made. Johnson County Crime Lab (JCCL) was one of the labs selected to do alpha testing on the kit and is currently using it for casework. The changes included a different incubation temperature (70°C vs. 56°C) and a plastic change in the plungers. These changes caused an overall improvement in yield. JCCL's organic procedure was compared to the "old" Casework sample kit and the "new" DNA IQ™ Casework Pro kit for Maxwell 16®. The percent change and total yield will be compared and discussed.