

Genomic DNA Purification from Cigarette Butts and Buccal Swabs Using the DNA IQ™ System

ABSTRACT

We describe experiments examining the ability of the DNA IQTM System to isolate DNA from cigarette butts and cigarette butt paper in the presence and absence of a proteinase K digestion. In addition we describe a protocol for DNA isolation from buccal swabs in the absence of a centrifugation step. These methods simplify the pretreatment steps for these samples and demonstrate that centrifugation of the lysate following initial incubation may not be required for successful DNA isolation. Elimination of centrifugation steps make complete processing of these sample types on automated workstations possible.



Download Article

(487 KB)