

INTERNAL VALIDATION OF KITS FOR USE IN TYPIFICATION OF REFERENCE SAMPLES

Alejandra Torres, Leticia Repetto, Juan Martín Marques, Laboratorio GeniaGeo

With the objective of having a large number of markers to establish reference samples without the need to extract and quantify the DNA from the same, is that tests were performed based on different samples in different media and treated with different reagents. Subsequently, following the recommendations of SWGDAM proceeded to internal validation kits; AmpFISTR® Identifiler® Direct, AmpFISTR® NGM Select™ Express and PowerPlex® 16 System, using Whatman FTA™ blood CARD and COPAN NUCLEIC-Card™ and oral mucosa obtained with brushes which were treated with Prep'n-Go™ Buffer and Swab Solution™ prior to amplification.