


# Validation Certificate

## VersaPlex® 27PY System

Product: DC7020

Developmental validation experiments and concordance studies were performed by Promega Corporation in collaboration with external laboratories to validate the performance of the VersaPlex® 27PY System for forensic purposes according to the guidelines set forth by the DNA Advisory Board (Quality Assurance Standards for Forensic DNA Testing Laboratories, 2020) and the Scientific Working Group on DNA Analysis Methods (Validation Guidelines for DNA Analysis Methods, 2016). The results of these validation experiments confirmed that this system can successfully be used for human identification testing by forensic and paternity laboratories. Each laboratory using the VersaPlex® 27PY System should perform its own appropriate internal validation studies to establish interpretation criteria and demonstrate that the VersaPlex® 27PY System is appropriate and fit for its own human identification uses.

Issued: February 2021  
Document #: VCDC70XX0221  
Promega Corporation  
2800 Woods Hollow Road  
Madison, WI 53711-5399  
United States

  
Douglas R. Storts  
Head of Research