



Promega

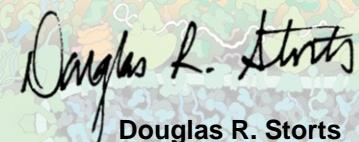
# Validation Certificate

## PowerPlex® Fusion System

Products: DC2402 and DC2408

Developmental validation experiments were performed by Promega Corporation in collaboration with external laboratories to validate the performance of the PowerPlex® Fusion System for forensic purposes according to the guidelines set forth by the DNA Advisory Board (Quality Assurance Standards for Forensic DNA Testing Laboratories, 2011) and the Scientific Working Group on DNA Analysis Methods (SWGDAM, Validation Guidelines for DNA Analysis Methods, 2012). The results of these validation experiments confirmed that this STR system can successfully be used for human identification testing by forensic and paternity laboratories. Each laboratory using the PowerPlex® Fusion System should perform its own appropriate internal validation studies to establish interpretation criteria and demonstrate that the PowerPlex® Fusion System is appropriate and fit for its own human identification uses.

Issued: November 2014  
Document #: VCDC24XX1114  
Promega Corporation  
2800 Woods Hollow Road  
Madison, WI 53711-5399  
United States



Douglas R. Storts  
Head of Research