

VALIDATION OF PLEXOR HY SYSTEM ON BIOMEK 3000

A. Abati

University of Liege, Forensic Pathology, Liege, Belgium

We'll present the validation performed for the validation of Plexor HY System on our Biomek 3000.

The validation work consist with establishing the best protocol to obtain the best Autosomal and Y Standard curves as possible.

After we have established the liquid handling best protocol, we verify the repeatability and repetitivity of the results, before using the Standard curve and after samples. We also compare the results with Quantifiler results on forensic samples.

We'll give all the parameters use for programming the Biomek 3000 and we'll show the results obtained before we decided to use Plexor HY System on Biomek 3000 for our forensic samples.