



European Forensic DNA Working Group Meeting

8th - 10th of November 2016, Florence, Italy

Agenda

Tuesday, 8th of November 2016

19.00 **Welcome Reception for networking at the hotel, followed by dinner**

Wednesday, 9th of November 2016

- 9.00 – 9.10 Welcome
Stephen Lupton, Promega Corporation, Paris, France
- 9.10 – 10.30 Session: New kits — new procedures**
- 9.10 – 9.30 Validation of the PowerQuant® system
Dave Moore, LGC, Teddington, UK
- 9.30 – 9.50 Prediction of autosomal STR typing success in ancient and Second World War bone samples
Irena Zupanic Pajnic, Institute of Legal Medicine, Ljubljana, Slovenia
- 9.50 – 10.10 Normalization and Direct Amplification of Casework Samples
Massimiliano Stabile, Forensic Scientist, Parma, Italy
- 10.10 – 10.30 The crash of FlyDubai Flight FZ981: a DNA approach to identification of vast body fragments
Pavel Ivanov, Federal Center for Forensic Medical Expertise of Ministry of Health, Moscow, Russia
- 10.30 – 11.00 Coffee break**
- 11.00 – 12.30 Session: NGS**
- 11.00 – 11.20 Next Generation Sequencing in Forensics: Achievements and Perspectives
Walther Parson, Institute of Legal Medicine, Innsbruck, Austria
- 11.20 – 11.40 SNPs, STRs and SNP-STRs - what genetic data can we get from Forensic MPS
Chris Philips, University of Santiago de Compostella, Spain
- 11.40 – 12:00 The Impact of Massively Parallel Sequencing on Genomics and a Vision for Forensic Genomics
Prof. Peter de Knijff, Human Genetics LUMC, Leiden, The Netherlands
- 12.00 – 12.30 Exploring sequence variation in forensic markers using next-generation sequencing and the PowerSeq® Auto/Mito/Y System
Tunde Huszar, University of Leicester, Alec Jeffreys Forensic Genomics Unit, Leicester, UK

12.30 – 14.30	Lunch
14.30 – 16.00	Breakout Session – Spectrum CE Instrument (Instrument, Software, Chemistry, Service)
16.00 – 17.00	Coffee break with Poster Session
18.30	Meet at hotel reception for evening event or at 19.00 in the restaurant “Il teatro del Sale”

Thursday, 10th of November 2016

9.00 – 9.20	Mass Disaster: DNA - a powerful tool for identification <i>Sophie Liénart, IFEG, Nantes, France</i>
9.20 – 9.40	PowerPlex® Y23 casework pilot study <i>Jim Thomson, LGC, Teddington, UK</i>
9.40 – 10.00	Y-STRs for paternity and clinical applications <i>Caroline Turner, Lancet, Johannesburg, South Africa</i>
10.00 – 10.20	Findings from the UK YHRD expansion project <i>Denise Syndercombe-Court, King's College London (KCL), London, UK</i>
10.20 – 10.40	Can some F.A.T. make your validation smoother? <i>Didier Hatsch, BioSilicium, Saint-Beauzire, France</i>
10.40 – 11.30	Coffee break and time for check out
11.30 – 12.00	Troubleshooting, Tips & Tricks <i>Rita Weispfenning, Promega Corporation, Stuttgart, Germany</i>
12.00 – 12.45	What's New from Promega? <i>Ann MacPhetridge, Promega Corporation, Madison, USA</i>
12.45 – 13.00	Discussion and Wrap-up
13.00 – 14.00	Lunch
14.00	End of meeting





European Forensic DNA Working Group Meeting

8th - 10th of November 2016, Florence, Italy

Poster

- 1) **NGS bacterial 16S RNA analysis in forensic soil identification**
Jakub Czarny, Institute of Forensic Genetics, Bydgoszcz, Poland
- 2) **Forensic SNP genotyping using array-based DNA Imaging Technology**
Jakub Czarny, Institute of Forensic Genetics, Bydgoszcz, Poland
- 3) **Application of an NGS-based age prediction model**
Anastasia Aliferi, King's College London (KCL), London, UK
- 4) **Development of PowerSeq™ Systems for Forensic Identification Using Massively Parallel Sequencing**
Spencer Hermanson, Margaret Ewing, Robert S. McLaren, and Douglas R. Storts
Promega Corporation, 2800 Woods Hollow Road, Madison, WI 53711, U.S.A.
- 5) **Y-Screening and Direct Amplification of Sexual Assault Samples**
Jeanne Bourdeau-Heller¹, Robert S. McLaren¹, Amy B. McGuckian², Jerome G. Remm², Julie Conover Sikorsky² and Douglas R. Storts¹.
¹Promega Corporation, ²Palm Beach County Sheriff's Office
- 6) **Normalization and Direct Amplification of Casework Samples**
Jeanne Bourdeau-Heller¹, Robert S. McLaren¹, Massimiliano Stabile², Nicola Staiti², Stella Eugenia Cirillo³, Lago Giampietro², and Douglas R. Storts¹.
¹Promega Corporation, ²Sez. Biologia, Parma, Italy, ³Università di Milano Bicocca
- 7) **Meeting the New Forensic ISO Standard: ISO 18385:2016**
Kristina Pearson and Susan Wigdal, Promega Corporation
- 8) **Comparing PowerPlex kit performance on Eppendorf vs 9700 cyclers**
PCR OPTIMIZATION OF STR SYSTEMS ON MULTIPLE THERMAL CYCLERS
Doug Wieczorek, Ph.D., Robert McLaren, Ph.D., Lotte Downey, and Rohaizah James, Ph.D., Promega Corporation
- 9) **Development Studies carried out for the processing of Crime Stain Samples on the IntegenX RapidHIT®200 platform**
Jenny Pagram, Lesley Ives, Key Forensic Services Ltd, Wymondham, UK
- 10) **Protocol Development for the IntegenX® RapidHIT® ID for Profiling Subject Samples**
D. Shackleton, N. Gray, D. King, S. Salceda, J. Buscaino, Lesley Ives, Key Forensic Services Ltd, Wymondham, UK
- 11) **NGS approach for the analysis of highly degraded bone samples**
Elena Pilli¹, Stefania Vai¹, Vincenzo Verdoliva², Marco Gigliucci², David Caramelli¹, Andrea Bert²,
¹Molecular Anthropology and Forensic Unit, University of Florence, Florence, Italy
²Sezione di Biologia, Rome, Italy

- 12) **PowerPlex® Fusion 6C contribution to probabilistic deconvolution of complex forensic samples**
Filippo Barni¹, Vincenzo Verdoliva¹, Andrea Berti¹, ¹Sezione di Biologia, Rome, Italy
- 13) **Evaluation of PowerQuant® System Accuracy in Complex Forensic Caseworks**
Simona Severini¹, Federica Tommolini², Gabriele Margiotta¹, Eugenia Carnevali¹
¹Department of Surgical and Biomedical Sciences, Section of Legal Medicine, Univeristy of Perugia, Perugia, Italy
²Department of Medical Biotechnologies, University of Siena, Siena, Italy; Department of Surgical and Biomedical Sciences, Section of Legal Medicine, Univeristy of Perugia, Perugia, Italy
- 14) **Irisplex System: Preliminary Findings of a Populationistic Study on an Umbrian Sample**
Federica Tommolini¹, Simona Severini², Gabriele Margiotta², Eugenia Carnevali²
¹Department of Medical Biotechnologies, University of Siena, Siena, Italy; Department of Surgical and Biomedical Sciences, Section of Legal Medicine, University of Perugia, Perugia, Italy
²Department of Surgical and Biomedical Sciences, Section of Legal Medicine, University of Perugia, Perugia, Italy
- 15) **Investigation into the sequence structure of 23 autosomal STRs and 23 Y-STRs using massively parallel sequencing**
So Yeun Kwon^{1,2}, Eun Hye Kim¹, Eun Young Lee¹, Hwan Young Lee¹, Kyoung-Jin Shin^{1,2}
¹Department of Forensic Medicine, Yonsei University College of Medicine, Seoul, Korea
²Brain Korea 21 PLUS Project for Medical Science, Yonsei University, Seoul, Korea
- 16) **DNA Identification of the victims from Italian aircraft crash near Skopje**
Zlatko Jakovski, Biljana Janeska, Natasa Davceva, Renata Jankova-Ajanovska, Ljupco Cakar, Natasa Bitoljanu, Violeta Belokaposka, Ana Ivceva, Sase Risteski, Hilda Jovanovic, Marija Bujaroska, Klimentina Trajkova, Aleksej Duma
Institute of Forensic Medicine, Criminology and Medical Deontology, School of medicine, Univ. Sn't Cyril and Methody, Skopje, Republic of Macedonia
- 17) **Discovery of skeletal remains in Flumineddu's cave in Sardinian 'Supramonte' (Italy): genetics's contribution to multidisciplinary reconstructive identification.**
Nutini A.L.¹, Fiaccabrino S.¹, Ricci U.¹, Carboni I.¹, Iozzi S.¹, Pinchi V.³, Focardi M.³, Norelli GA.³ Piredda V.², Mazzeo E.²
¹ SOD Genetic Diagnostic Azienda Ospedaliera-Universitaria Careggi of Florence, Italy
² Department of Biomedical Sciences, Legal Medicine, University of Sassari, Italy
³ Department of Health Sciences, Forensic Sciences Section, University of Florence, Italy

