



Mr. Michael Spiegel
Deputy Director MATF, Monash University (Australia)

Biography

Michael has worked in the Life Sciences arena for 13 years, including eight years at the European Molecular Biology Laboratory. He has worked in the laboratory of the EMBL Director General, Fotis Kafatos in Heidelberg Germany and in the antibody core facility with Alan Sawyer. Michael has managed facilities studying immune system and antibody production, and spent a number of years in the EMBL animal facilities unit. He has extensive experience managing large teams of diverse individuals, has a proven track record of tackling technologically difficult projects taking them from idea to implementation. He is currently deputy director of the recently established Monash Antibody Technology Facility at Monash University.