

EXPLORING ZOOONOTIC PATHOGENS IN AUSTRALIAN LIVESTOCK.

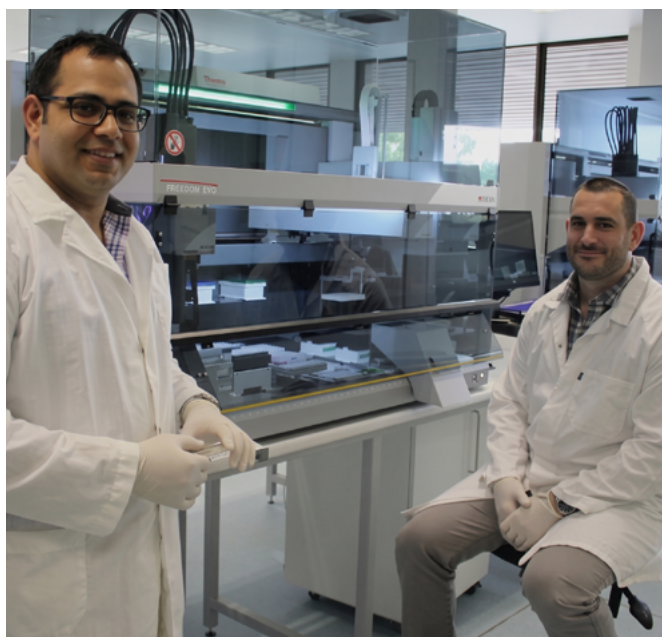
Tecan Journal article.



Antimicrobial resistance (AMR) is a growing issue preventing the effective treatment of numerous parasitic diseases in both humans and animals. The lack of effective antimicrobials for the treatment of a growing number of diseases represents a serious threat to global public health, making it imperative that governments and societies around the world take action to tackle this issue. Scientists at the University of Murdoch in Australia are studying antimicrobial resistance in livestock – which can potentially spread to humans – using a statistical approach of surveillance to monitor the situation more closely.

Antimicrobial resistance is causing a growing number of standard antibiotic, antifungal, antiviral and antimalarial treatments to become ineffective, leading to increased prevalence and spread of many infectious diseases. This issue is not isolated to human medicine and is a growing concern for veterinary medicine. It is particularly problematic in livestock, due to a combination of relatively concentrated animal populations and the rise of drug-resistant zoonotic pathogens that can spread between animals and humans.

Dr Sam Abraham, a microbiologist at Murdoch University in Perth, Australia, is investigating AMR at the interface between human and animal health, primarily by studying zoonotic pathogens that are present in Australian livestock. Sam explained: “Antimicrobial resistant bacteria that emerge and spread in livestock and other animals contribute to the global burden of AMR. As a result, there is a significant public health concern over the transfer of such bacteria to humans via food and the environment and the rapid transmission of the genes responsible for AMR into other human pathogens.”



Tecan's automated solutions and library preparation kits are helping Sam Abraham (left) and his colleagues to study antimicrobial resistance in livestock.

“Our lab focuses on determining the minimum inhibitory concentration (MIC) of drugs for bacteria found in fecal samples from chickens, pigs, sheep and cattle – such as *E. coli*, *Salmonella*, *Staphylococcus* and *Streptococcus* – and characterizing the genomic mechanisms of drug resistance in these bacteria. This project is primarily funded by the Australian government, but we also have global collaborators. The aim of this work is to create a snapshot database of populations of enteric bacteria and, once optimized, use it to monitor levels of antibiotic resistance in livestock all over the world.¹ We use a range of techniques, including classical antibiotic susceptibility testing (AST), RT-PCR, NGS, MALDI-TOF and traditional bacteriology methodologies, such as gram staining.”

“This type of surveillance project requires large-scale sampling of indicator organisms to identify AMR and monitor whether these resistances are being detected more frequently,” Sam continued. “Unfortunately, most countries only use low numbers of animal samples for surveillance, which is unlikely to highlight emerging issues until it is too late. This is because conventional, human-centric surveillance methods are hamstrung by high processing costs and slow turnaround times, making them incompatible with the high volumes of sampling required to accurately depict a population's AMR status. We therefore need to use high throughput automation and genomics solutions with rapid turnaround times, enabling the analysis of large data sets to help identify emerging problems quickly and respond to those problems rapidly.”

“I was initially introduced to Tecan when I was a postdoc setting up the first national network on AMR surveillance of Australian animals. I saw one of its trade displays for microbiology and chatted to the very knowledgeable team, who helped us to select a solution that would best suit our AMR screening requirements. In 2016, we were able to purchase our Tecan systems using grants from the Department of Agriculture and Australian Pork Limited. The Tecan

team here in Australia was pivotal to this project, partnering with us on the grant application, and working closely with us to design an automation set-up that would meet our needs. We currently have two Freedom EVO® 150 workstations, one of which has a number of SciRobotics devices integrated onto it to create our Robotic Antimicrobial Susceptibility Platform (RASP).²

“Our workflow starts with a homogenized sample, which is transferred to the Freedom EVO – where it could potentially undergo various dilutions depending on the sample – and automatically plated out using the PetriPlater™. The plates are stacked and stored in a PetriSel™ and all of our tubes and plates are barcoded so that the samples can be automatically tracked. This makes the whole system very efficient and troubleshooting is much easier if there are ever any problems. We plate on chromogenic agar, which allows our PickoloMI™ robotic colony picker – programmed with assistance from Tecan – to easily identify and differentiate specific micro-organisms. These colonies are then picked and grown – either for a few hours or overnight – in a 96-well plate.”

“The number of errors that we generate is also hugely minimized – the beauty of automation – and this results in excellent quality data.”

“From here, there are different routes of analysis depending on the sample and what we are looking for. Some we send to MALDI-TOF mass spectrometry analysis for bacterial identification or we perform AST in house to determine the AMR profile. We have also integrated next generation sequencing with our microbiology workflow, using a Freedom EVO 150 to perform automated library preparation with Tecan’s Celero™ EZ DNA-Seq library preparation kit. This allows us to rapidly perform genotyping or whole genome sequencing on bacteria that have

demonstrated AMR. Genotyping gives us the full picture about the resistant pathogen: is it human-, livestock- or even gull-derived? Is it only a problem for animal health, human health or both? We can answer all these questions with genomics. All the data generated is then sent to and stored in our national database.”

“Using automation, we can process 5,000 to 10,000 bacterial isolates a week – around 150 plates a day – compared to a couple of hundred isolates manually. This means that we only need one researcher working on this at a time and, as our throughput has increased so significantly, others can spend time on different tasks, which is an incredibly effective time saver. The number of errors that we generate is also hugely minimized – the beauty of automation – and this results in excellent quality data, which is fantastic. We are very happy with the instruments,” Sam concluded.

Literature

- 1) Laird TJ, et al. J Antimicrob Chemother, 2021. doi: 10.1093/jac/dkab403. Online ahead of print.
- 2) Truswell A, et al. J Antimicrob Chemother, 2021, 76, 1800-1807. doi: 10.1093/jac/dkab107.

Additional resources

To explore Tecan’s veterinary solutions, visit:
www.tecan.com/veterinary

To find out more about Tecan and SciRobotics’ colony picking and plating solutions, visit:
www.tecan.com/colonypicking

To learn more about the Antimicrobial Resistance and Infectious Diseases Laboratory (AMRID) at Murdoch University, go to <https://amrid.com.au/>



For research use only. Not for use in diagnostic procedures.

.....
Australia +61 3 9647 4100 **Austria** +43 62 46 89 330 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170
India +91 12 4494 2050 **Italy** +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Nordic** +46 8 750 39 40 **Singapore** +65 644 41 886
South Korea +82 2 818 3301 **Spain** +34 93 595 25 31 **Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +41 44 922 81 11
.....

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication, however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. All mentioned trademarks are protected by law. In general, the trademarks and designs referenced herein are trademarks, or registered trademarks, of Tecan Group Ltd., Männedorf, Switzerland. A complete list may be found at <http://www.tecan.com/trademarks>. Product names and company names that are not contained in the list but are noted herein may be the trademarks of their respective owners. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative.

Tecan is in major countries a registered trademark of Tecan Group Ltd., Männedorf, Switzerland.

© 2026 Tecan Trading AG, Switzerland, all rights reserved.

www.tecan.com