



2016 Travel Award Application (cont.)

Please describe the role MEA data plays in your poster:	
Application Area	
Cell type	
Plate type	

Briefly describe your experimental design:

--

Briefly describe your MEA results:

--

How did MEA data influence your conclusions?

--

Email completed form to: marketing@axionbio.com