



Travel Award Application

Email completed form to: marketing@axionbio.com

Contact information:	
Name	
Organization	
Department/Lab	
Address	
Street	
City	
State/Province	
Postal Code	
Country	
Phone	
Email	

Tell us about the conference you'd like to attend:	
Conference	
Date	
Location	

Abstract (please attach a copy of your abstract acceptance from the conference organizers):

--

Please provide more information about your MEA experiments on Page 2.



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