

# Enter to win a Fitbit!

Visit both InSphero (BOOTH #113) and ALPCO (BOOTH #1649) for a chance to win a Fitbit.

- 1 Complete the contact information section of the entry form.
- 2 Ask a staff member in each booth to stamp your entry form.
- 3 When you have both stamps, detach the entry form at the perforation.
- 4 Place it in the entry box in either booth for your opportunity to win!

*One entry per person. The winner will be contacted via email and the Fitbit will be mailed to the address provided on the entry form.*

## CONTACT INFORMATION

Name

Company

Address

City, State, Zip

Phone

Email

## INSPHERO STAMP (BOOTH #113)

## ALPCO STAMP (BOOTH #1649)

Enter to win  
a Fitbit!



*See other side for details.*

Join us as Dr. Burcak Yesildag, InSphero AG, presents:

**POSTER 2202-P**

Using Uniform Reaggregated Pancreatic Islets in a Microfluidic Perifusion System Enables Studying Insulin Release Dynamics at Single-Islet Level

**GENERAL POSTER SESSION**

Sunday, June 11, 12:00 – 1:00 pm

**MODERATED POSTER DISCUSSION**

Monday, June 12, 12:00 – 1:00 pm

Booth #113



Booth #1649

