

Cancer and Metabolism 2018

Cambridge, UK
June 25 - 27, 2018

Organizer:

- Christian Frezza (MRC Cancer Unit, UK)
- Eyal Gottlieb (Israel Institute of Technology, Israel)
- Karen Vousden (The Francis Crick Institute, UK)

Speakers:

- Heather Christofk (UCLA, US)
- Ralph DeBerardinis (UT Southwestern, US)
- Aimee Edinger (UC Irvine, US)
- Martin Eilers (Würzburg Universität, Germany)
- Ayelet Erez (Weizmann Institute, Israel)
- Sarah-Maria Fendt (VIB, Belgium)
- Matt Vander Heiden (MIT, US)
- Oliver Maddocks (Beatson Institute, UK)
- Brendan Manning (Harvard University, US)
- Josh Rabinowitz (Princeton University, US)
- Alexei Vazquez (Beatson Institute, UK)
- Kathryn Wellen (University of Pennsylvania, US)

Talk and poster abstracts invited

- More information available on the meeting website