



European Laboratory  
Research & Innovation Group



Miltenyi Biotec



**TOCRIS**  
a biotechnique brand



ELRIG Networking Event in Association with AstraZeneca

**“PROTACs – An exciting therapeutic modality”**

6<sup>th</sup> December 2017, 14.00 – 19.00

**Alderley Park Conference Centre**

Alderley Park, Nether Alderley, Macclesfield, Cheshire, SK10 4TG

Proteolysis targeting chimeras (PROTACs) offer a novel alternative to traditional small molecules. They inhibit proteins by binding to their functional regions and by recruiting E3 ligases to target the protein for degradation through the cells natural ubiquitin/proteasome pathway. This event will focus on the recent developments in the field, provide an opportunity to learn about the types of protein complexes that are being considered as PROTAC targets and provide an introduction to those new to the field who are looking to start engaging with PROTACs within their own research. Informal presentations will be given by guest speakers followed by a networking session over food and drinks.

**Confirmed speakers for this event are:**

**Shaun Cowley** (Chair of Molecular Biology, University of Leicester)

**Thomas Hayhow** (Medicinal Chemist, AstraZeneca)

**John Spencer** (Professor of Bio-organic Chemistry, University of Sussex)

**Jason Brown** (Scientific Director, Ubiquigent)

**Why attend?**



**Communicate**

Expand your local network and share research with like-minded individuals



**Collaborate**

Explore collaboration opportunities between academia and the pharma/biotech industry



**Innovate**

Accelerate the pace of innovation through early development partnership

ELRIG networking events are FREE-TO-ATTEND.  
For further information or to register, contacting [booking@elrig.org](mailto:booking@elrig.org)  
Follow us on Twitter @ELRIG\_UK for regular updates

Coming soon - Look out for the dates for our 2018 networking events on the ELRIG website:  
[www.elrig.org/networking](http://www.elrig.org/networking)