

# 7<sup>th</sup> NEB Meeting on DNA Restriction & Modification

Aug. 24–29, 2015 | Univ. of Gdansk, Poland



UNIVERSITY OF GDANSK

The Seventh New England Biolabs Meeting on DNA Restriction and Modification will take place at the University of Gdansk, Poland, August 24-29, 2015. The agenda includes morning and afternoon/evening sessions; a continuous poster session; and free time for discussions. Sessions will be held at the Intercollegiate Faculty of Biotechnology in the historic part of the city near the old town center. The meeting web site accepts registrations, abstracts, and payments by credit card: <http://nebmeetinggdansk2015.pl>

Oral sessions will comprise 30-minute talks from invited speakers and shorter talks from attendees based on their abstracts. The first talk of each session will be longer and will also provide a general overview of field. Thanks to generous sponsorship, the cost of the meeting has been kept to a minimum. Scholarships are available for young scientists, and financial help for established scientists who cannot otherwise attend.

## Session Topics:

- Protein structures and interactions
- Genetics and biology of restriction, modification, and accessory enzymes
- Mechanisms of catalysis and DNA sequence-recognition
- Protein engineering and emerging technologies
- Gene targeting: zinc fingers, meganucleases, TALENs, CRISPR/Cas 9
- Epigenetics; bacterial methylomes; alternative DNA-modifications

## Confirmed Speakers:

Aneel Aggarwal	Mt. Sinai School of Medicine, USA
Jens Boch	Martin Luther University, Halle, Germany
Matthias Bochtler	Intl. Inst. Molecular & Cell Biology, Warsaw, Poland
Robert Blumenthal	University of Toledo College of Medicine, USA
Janusz Bujnicki	Intl. Inst. Molecular & Cell Biology, Warsaw, Poland
Toni Cathomen	Universitätsklinikum Freiburg, Germany
Srinivasan Chandrasegaran	Johns Hopkins School of Public Health, USA
Xiaodong Cheng	Emory University School of Medicine, USA
David Dryden	Edinburgh University, UK
Yuval Ebenstein	Tel Aviv University, Israel
David Edgell	University Western Ontario, Canada
Elizaveta Gromova	Moscow State University, Russia
Steve Halford	University of Bristol, UK
Nancy Horton	University of Arizona, USA
Petra Imhof	Freie Universität Berlin, Germany

Keiko Kita	Kyoto University, Japan
Ben Kleinstiver	Harvard Medical School, USA
Saulius Klimasauskas	Inst. of Biotechnology Vilnius University, Lithuania
Ichizo Kobayashi	University of Tokyo, Japan
Ite Laird-Offringa	University of Southern California, USA
Heinrich Leonhardt	Ludwig Maximilians University, Germany
Valakunja Nagaraja	Indian Institute of Science, Bangalore
Robert Neely	University of Birmingham, UK
Sriharsa Pradhan	New England Biolabs, Inc., USA
Desirazu Rao	Indian Institute of Science, Bangalore
Norbert Reich	U.C. Santa Barbara, USA
Nigel Richards	Indiana University-Perdue University, USA
Rich Roberts	New England Biolabs, Inc., USA
Virginijus Siksnys	Inst. of Biotechnology Vilnius University, Lithuania
Barry Stoddard	Fred Hutchinson Cancer Res. Center, USA
Elmar Weinhold	RWTH Aachen University, Germany
Guo-Liang Xu	Chinese Academy of Sciences, Shanghai

### Deadlines:

July 10:	Last day for regular registration. Later registrations incur a late fee
July 17:	Last day to apply for a young scientist scholarship.
July 24:	Abstracts sent after July 24 <sup>th</sup> are unlikely to be chosen for short talks.
July 31:	Abstracts sent after July 31 <sup>st</sup> are unlikely to appear in the Abstract Book.
Aug 24:	Last-minute, walk-in registration if full-capacity is not already reached.

### Registration (includes conference, lunches, dinners, tour and banquet):

	Academic/Commercial	Student/Post-Doc
Regular rate:	350 Euro	275 Euro
After July 10th:	400 Euro	300 Euro

### Accommodation:

Dormitory (5 nights):	100 Euro (double)	175 Euro (single)
Local Hotels: ~75-125 Euro per night. Please make your own arrangements		

**Organizers:** Iwona Mruk (Poland): [contact@nebmeetinggdansk2015.pl](mailto:contact@nebmeetinggdansk2015.pl)  
 Geoff Wilson (USA): [wilson@neb.com](mailto:wilson@neb.com)  
 Rick Morgan (USA): [morgan@neb.com](mailto:morgan@neb.com)



**Previous NEB meetings:** 1988: **Gloucester**, USA  
 1990: **Berlin**, Germany  
 1994: **Vilnius**, Lithuania  
 1997: **Igls**, Austria  
 2004: **Bristol**, England  
 2010: **Bremen**, Germany