

PROGRAM ER & REDOX CLUB MEETING 15-17 APRIL 2015 VENICE

APRIL 15th

14:00	14:20	WELCOME & OPENING ADDRESS	
14:20	15:40	SESSION 1: ENTER THE ER <i>chair: Ester Zito</i>	
14:20	14:40	Tsilil Ast (M. Schuldiner)	Forget the plunger - Unclogging the translocon
14:40	15:00	Stefan Schorr	Co-chaperone specificity in gating of the polypeptide conducting channel in the membrane of the human ER
15:00	15:20	Matteo Fossati (N. Borgese)	A positive signal restrains rab6-dependent recycling of membrane proteins in the early secretory pathway
15:20	15:40	Luca Scorrano	Why do we have the ER? A mitochondrial answer
15:40	16:00	Coffee break	
16:00	17:20	SESSION 2: ER CHAPERONES <i>chair: Roberto Sitia</i>	
16:00	16:20	Matthias Feige	The sequence-based quality control code of the Hsp70 chaperone system in the endoplasmic reticulum
16:20	16:40	Mathias Rosam (J. Buchner)	BAP, an atypical NEF for BiP
16:40	17:00	Naomi Lodder (I. Braakman)	Characterising the function of pERp1
17:00	17:20	Alessandro Bosetti (Promega)	Interrogating protein dynamics within living cells using protein-based reporters derived from NanoLuc Luciferase
17:20	17:45	Aperitivo	
17:45	18:30	1st POSTER ADVERTS SESSION <i>chair: Eelco van Anken</i>	
18:30	20:00	Dinner	
20:00	22:00	1st POSTER SESSION	

APRIL 16th

7:30	9:00	Breakfast	
9:00	10:20	SESSION 3: PDI FAMILY <i>chair: Neil Bulleid</i>	
9:00	9:20	William Remelli (A. Ceriotti)	Characterization of recombinant AtPDL5-1, the smallest putative PDI-like protein of Arabidopsis thaliana
9:20	9:40	Helena Safavi-Hemami (L. Ellgaard)	Specialized venom gland protein disulfide isomerases facilitate the folding of cone snail toxins
9:40	10:00	Giorgia Brambilla Pisoni (M. Molinari)	Division of labor between oxidoreductases: TMX1 preferentially acts upon membrane-bound polypeptides
10:00	10:20	Tiziana Anelli (R. Sitia)	Proteostasis and redoxstasis in the secretory pathway: tales of tails from ERp44 and immunoglobulins
10:20	10:40	Coffee break	
10:40	12:00	SESSION 4: DISULFIDE BOND FORMATION <i>chair: Aldo Ceriotti</i>	
10:40	11:00	Lei Wang (C. Wang)	Interplay between Ero1 and PDI on oxidative protein folding in the endoplasmic reticulum
11:00	11:20	Adam Benham	Exploring the effect of the redox environment on Ero1 & PDI
11:20	11:40	Philip Robinson (N. Bulleid)	Protein structure and disulphide formation following nascent chain exposure to the ER lumen
11:40	12:00	Lloyd Ruddock	Applications of disulfide bond formation in the cytoplasm
12:30	14:00	Lunch	
14:20	15:40	SESSION 5: UNFOLDED PROTEIN RESPONSE <i>chair: Luca Rampoldi</i>	
14:20	14:40	Florentina Pena	The endoplasmic reticulum under hypoxic stress
14:40	15:00	Tomás Aragón	A new mechanism for XBP1 mRNA transport to ER stress signaling centers
15:00	15:20	Andrea Orsi (E. van Anken)	ATF6 and ERAD are key for transitioning from acute to chronic ER stress responses
15:20	15:40	David Ron	Further molecular insight into the working of the Integrated Stress Response, ISR
15:40	16:00	Coffee break	
16:00	17:20	SESSION 6: ER-STASIS IN HEALTH & DISEASE <i>chair: David Ron</i>	
16:00	16:20	Carmine Settembre	Self eating during vertebrate development
16:20	16:40	Matteo Trudu (L. Rampoldi)	ER stress in models of uromodulin-associated kidney disease
16:40	17:00	Vera Volpi (M. D'Antonio)	The role of the ERAD pathway in the physiology and disease of peripheral myelination
17:00	17:20	Ester Zito	ER-localized selenoprotein SEP1, linked to skeletal muscle pathology, redox-regulates SERCA2 pump activity
17:20	17:45	Aperitivo	
17:45	18:30	2nd POSTER ADVERTS SESSION <i>chair: Johannes Buchner</i>	
18:30	20:00	Dinner	
20:00	22:00	2nd POSTER SESSION	
22:00	23:30	OPTIONAL: trip "Venice by night" to Piazza San Marco & Basilica	

APRIL 17th

7:30	9:00	Breakfast	
9:00	10:20	SESSION 7: ER-ASSOCIATED DEGRADATION <i>chair: Maurizio D'Antonio</i>	
9:00	9:20	Ryo Ushioda (K. Nagata)	ERdj5-mediated ER homeostatic mechanism
9:20	9:40	Kenji Inaba	Dual conformations of ERdj5 play a significant role in acceleration of the ER associated degradation
9:40	10:00	Federica Lari (J. Christianson)	UPRegulating ERAD: role of ER-membrane ubiquitin ligases in ER homeostasis
10:00	10:20	Anett Köhler (T. Sommer)	The role of the mannosidase Htm1p in glycoprotein ERAD
10:20	10:40	Coffee break	
10:40	11:20	SESSION 8: ER-ASSOCIATED DEGRADATION <i>chair: Maurizio Molinari</i>	
10:40	11:00	Stephen High	Small glutamine-rich tetratricopeptide repeat-containing protein alpha (SGTA), saint or sinner?
11:00	11:20	Marius K. Lemberg	Sampling the membrane: function of intramembrane proteases in ER protein homeostasis
11:20	11:40	Ludger Johannes	Identification of cellular targets of Retro-2 that are required for retrograde toxin trafficking
11:40	12:15	CLOSING REMARKS: <i>chair: Ineke Braakman</i>	
		Awards, Evaluation & Election new organizers	
15:00	18:00	OPTIONAL: "boat trip through the lagoon"	

POSTER PRESENTATIONS

APRIL 15th

Anush Bakunts (E. van Anken)	<i>A double-edged sword: use of the APEX-tag for assessing changes in the morphology & proteome of the ER</i>
Nina Bergbold (M.K. Lemberg)	<i>The ERAD rhomboid protease RHBDL4 interacts with the ERAD machinery to modulate IP3R abundance</i>
Stefano Bestetti (R. Sitia)	<i>Regulation of H2O2 membrane transport during cell stress</i>
Anna Paola Casazza (A. Ceriotti)	<i>Preliminary characterization of AtPDI5-1</i>
Fiona Chalmers (N. Bulleid)	<i>Increasing the capacity of the ER to fold antibodies by manipulation of stress response pathways</i>
Anna Chatsivili (I. Braakman)	<i>PDI: how important for folding is the most abundant oxidoreductase of the ER?</i>
Ruming Chen (D. Ron)	<i>G-actin provides substrate-specificity to eukaryotic initiation factor 2α holophosphatases</i>
Mari Chiritoiu (S.M. Petrescu)	<i>The intrinsically disordered domain of EDEM1 is key for association with its clients</i>
Alberto Danieli (E. van Anken)	<i>A novel approach for correlative light and electron microscopy: the dark side of the μ</i>
Emma Fenech (J. Christianson)	<i>Constructing the ER E3 ubiquitin ligase interaction network</i>
Sefi Geva (M. Schuldiner)	<i>A systematic screen uncovers a novel cargo receptor</i>
Christine John (J. Buchner)	<i>Analysis of the interactions of variable antibody domains</i>
Shingo Kanemura (K. Inaba)	<i>A newly identified disulfide bond in regulation of human Ero1α</i>
Raffaella Magnoni (L. Ellgaard)	<i>FICD is a new unfolded protein response target involved in BiP AMPylation</i>
Cristina Scapin (M. D'Antonio)	<i>Translational attenuation via eIF2α-phosphorylation is protective in a POS63del-CMT1B neuropathic mouse model</i>

APRIL 16th

Marianna Marino (E. Zito)	<i>Functional interplay between SEPN1 and ascorbic acid</i>
Kitty McCaffrey (I. Braakman)	<i>The ER and Inflammation</i>
Daniel Mesquita da Fonseca (I. Braakman)	<i>CFTR NBD1 folding requires Hsp90</i>
Déborah Naón (L. Scorrano)	<i>Mitofusin 2 is alternatively spliced in ER specific variants that control ER-mitochondria tethering and ER shape</i>
Masaki Okumura (K. Inaba)	<i>Real-time monitoring of PDI-catalyzed oxidative protein folding by high-speed atomic force microscopy</i>
Džana Pašalić (J. Buchner)	<i>The tail piece dictates the oligomerization of IgM in vitro and in vivo</i>
Riccardo Ronzoni (A. Fra)	<i>Aberrant disulphide interactions contribute to the secretory deficiency of Alpha 1-Antitrypsin mutants</i>
Sara Sannino (R. Sitia)	<i>Dual regulation of ERp44 in the early secretory pathway</i>
Céline Schaeffer (L. Rampoldi)	<i>Identification and characterisation of cellular pathways activated or dysregulated by mutant uromodulin expression</i>
Emilie Schrepfer (L. Scorrano)	<i>Dissecting Mitofusin 2 interactors at the ER-mitochondria interface</i>
Chloe Stoylo (N. Bulleid)	<i>Antibody assembly and secretion in CHO cells</i>
Christina Stutzer (J. Buchner)	<i>Characterization of the surrogate light chain</i>
Milena Vitale (E. van Anken)	<i>Dynamics of endoplasmic reticulum stress responses to sudden and persistent proteostatic challenges</i>
Laura Tadè (E. van Anken)	<i>Homeostatic readjustment of the endoplasmic reticulum under chronic stress</i>
Uri Weill (M. Schuldiner)	<i>Swap-N-Tag: a new tool for systematic characterization of yeast proteins</i>
Cristina Scapin (ERSD)	<i>Endoplasmic Reticulum Stress in Diseases, a new journal focusing on the connection between ER stress & disease</i>

PARTICIPATING GROUPS

Tomás Aragón	<i>Centro de Investigación Médica Aplicada</i>	Pamplona	Spain
Adam Benham	<i>Durham University</i>	Durham	England
Nica Borgese	<i>Consiglio Nazionale delle Ricerche Milano</i>	Milan	Italy
Ineke Braakman	<i>Universiteit Utrecht</i>	Utrecht	Holland
Johannes Buchner	<i>Technische Universität München</i>	Munich	Germany
Neil Bulleid	<i>University of Glasgow</i>	Glasgow	Scotland
Aldo Ceriotti	<i>Consiglio Nazionale delle Ricerche Milano</i>	Milan	Italy
John Christianson	<i>Ludwig Cancer Research</i>	Oxford	England
Maurizio D'Antonio	<i>Istituto Scientifico San Raffaele</i>	Milan	Italy
Lars Ellgaard *	<i>Københavns Universitet</i>	Copenhagen	Denmark
Matthias Feige	<i>Technische Universität München</i>	Munich	Germany
Anna Fra	<i>Università degli studi di Brescia</i>	Brescia	Italy
Stephen High	<i>University of Manchester</i>	Manchester	England
Kenji Inaba	<i>Tohoku University</i>	Sendai	Japan
Ludger Johannes	<i>Institut Curie</i>	Paris	France
Marius K. Lemberg	<i>Zentrum für Molekulare Biologie</i>	Heidelberg	Germany
Maurizio Molinari	<i>Istituto di Ricerca in Biomedicina</i>	Bellinzona	Switzerland
Kazuhiro Nagata *	<i>Kyoto Sangyo University</i>	Kyoto	Japan
Florentina Pena	<i>Institutul de Biochimie al Academiei Române</i>	Bucharest	Romania
Stefana M. Petrescu *	<i>Institutul de Biochimie al Academiei Române</i>	Bucharest	Romania
Luca Rampoldi	<i>Istituto Scientifico San Raffaele</i>	Milan	Italy
David Ron	<i>University of Cambridge</i>	Cambridge	England
Anna Rubartelli	<i>Istituto Scientifico San Raffaele</i>	Milan	Italy
Lloyd Ruddock	<i>University of Oulu</i>	Oulu	Finland
Maya Schuldiner *	<i>Weizmann Institute</i>	Rehovot	Israel
Luca Scorrano	<i>Università di Padova</i>	Padua	Italy
Carmine Settembre	<i>Telethon Institute of Genetics and Medicine</i>	Pozzuoli	Italy
Roberto Sitia	<i>Istituto Scientifico San Raffaele</i>	Milan	Italy
Thomas Sommer *	<i>Max Delbrück Centrum</i>	Berlin	Germany
Eelco van Anken	<i>Istituto Scientifico San Raffaele</i>	Milan	Italy
Chih-chen Wang *	<i>Institute of Biophysics Chin. Acad. of Sciences</i>	Beijing	China
Richard Zimmermann	<i>Universitätsklinikum des Saarlandes</i>	Homburg	Germany

* Not personally present; representatives from their groups are