

Poster Index

- Poster 1 Evi Aerts (University of Antwerp, Belgium)
Identification of new missense variations in the SH2B1 gene in obese and lean individual
- Poster 2 Azizah Alanazi (King Faisal Specialist Hospital and Research Center, Saudi Arabia)
Inflammatory mediators in human adipocytes: Role of hypoxia
- Poster 3 Maria Alfaradhi (University of Cambridge, UK)
Oxidative stress and altered lipid homeostasis accompany fatty liver programmed by maternal obesity
- Poster 4 Neli Atanassova (Institute of Metabolic Sciences, UK)
KSR2 mutations cause obesity and insulin resistance by disrupting food intake, energy expenditure and substrate utilisation in humans
- Poster 5 Pernille Barkholt (Gubra ApS, Denmark)
Metabolic changes in rats following roux-en-y gastric bypass - peripheral and central modes of action
- Poster 6 Lindsey Berends (University of Cambridge, UK)
Early-life programming of central insulin sensitivity in adult male mice
- Poster 7 Luke Burke (University of Cambridge, UK)
5-HT obesity treatment anorectic efficacy is not compromised by ageing
- Poster 8 Francesca Cirulli (Istituto Superiore di Sanità, Italy)
The p66Shc -/- mouse as an animal model to study basic mechanisms regulating energy metabolism and its effects on healthspan
- Poster 9 Barbora Doslikova (University of Cambridge, UK)
Pro-opiomelanocortin neurons in the nucleus of the solitary tract are responsive to leptin and 5-HT
- Poster 10 Denise Fernandez-Twinn (University of Cambridge, UK)
Adipose tissue insulin signalling via IRS-1 is impaired in offspring of obese mice and is associated with sequestration by miR-126
- Poster 11 Katharina Gapp (Brain Research Institute Zurich, Switzerland)
Implication of sperm miRNAs in the transgenerational effects of early traumatic stress in mammals
- Poster 12 Clemente Garcia-Rizo (Fundacio Clinic per a la Recerca Biomedica, Spain)
Energy imbalance in cannabis users with a first episode of psychosis
- Poster 13 Michelle Goldsworthy (MRC Harwell, UK)
A recessive mouse mutagenesis ageing screen for metabolic and body composition phenotypes
- Poster 14 Safiya Hafeji (University of St Andrews, UK)
A pilot study investigating the influence of visceral obesity on outcomes following anterior resection complicated by anastomotic leak
- Poster 15 Nicholas Heeley (Institute of Metabolic Science, UK)
5-HT₆ receptor antagonist induced appetite suppression involves hypothalamic circuitry

- Poster 16 Andrew Holmes (University of Liverpool, UK)
RNAseq and histological analyses of hypothalami of lean and hypermetabolic Gnasxl knockout mice reveals reduced levels of the astrocyte marker Gfap
- Poster 17 Alison Hough (MRC, UK)
Functional analysis of the type 2 diabetes GWAS gene KLF14 using mouse models
- Poster 18 Camilla Ingvorsen (DTU-Systems Biology, Denmark)
Effect of gestation on obesity-induced hepatic and placental inflammation in mice
- Poster 19 Michelle Johnson (The Open University, UK)
Changes in appetite hormones and gut tissue architecture during the rat reproductive cycle, pregnancy and lactation
- Poster 20 Toshiya Kojima (University of Tokyo, Japan)
SUMOylation of Krüppel-like factor 5 controls food intake by suppressing AgRP expression on contact with FoxO1 in hypothalamic neurons
- Poster 21 Andreas Kolb (University of Aberdeen, UK)
Alpha casein deficiency attenuates growth in early postnatal life and extends life-span
- Poster 22 Sander Kooijman (Leiden University Medical Center, The Netherlands)
The central melanocortin system regulates VLDL-triglyceride clearance
- Poster 23 Jana Kracmerova (Charles University in Prague, Czech Republic)
Effect of acute increase of free fatty acids levels by intralipid infusion on inflammatory status of obese patients
- Poster 24 Jose Lopez-Pedrosa (Abbott Nutrition R&D, Spain)
Slow digesting carbohydrates during pregnancy alters genomic imprinting for enhancing muscle development in the offspring
- Poster 25 Marcella Ma (Institute of Metabolic Science, UK)
FTO interacts with RNA and its deficiency is associated with an increased burden of 3 methyluridine in RNA in vivo
- Poster 26 Adil Mardinoglu (Chalmers University of Technology, Sweden)
Global analysis of human adipocyte metabolism in response to obesity - a systems biology approach
- Poster 27 William Marsh (Metabolic Research Laboratories, UK)
Recurrent hypoglycaemia increases miRNA expression in the anterior cingulate cortex of brain
- Poster 28 Pablo B Martínez de Morentin (University of Santiago de Compostela, Spain)
Central actions of estrogens on energy balance
- Poster 29 Malgorzata Martin-Gronert (University of Cambridge, UK)
Global transcriptional profiling reveals 5-HT2AR as a hypothalamic target of developmental programming
- Poster 30 John Menzies (University of Edinburgh, UK)
Reciprocal interactions of stress and palatable food reward in prenatally stressed rats
- Poster 31 Myrte Merkestein (University of Oxford, UK)
FTO-induced changes in gene expression and RNA methylation in mice
- Poster 32 Lee Moir (MRC Harwell, UK)
An enu induced missense mutation in the transcription factor orthopedia homolog results in a novel mouse model of obesity

- Poster 33 Nicolas Nunn (University of Manchester, UK)
The thermogenic effects of leptin are mediated by hypothalamic PrRP neurones
- Poster 34 Dorota Olszewska-Slonina (Nicolaus Copernicus University in Torun, Poland)
Serum paraoxonase (PON155) polymorphism and PON1 activity in patients with different degree of obesity
- Poster 35 Edwin Parlevliet (Leiden University Medical Center, The Netherlands)
*The GLP-1 receptor agonist exendin-4 reduces non-alcoholic steatohepatitis in addition to atherosclerosis development in APOE*3.Leiden.CETP mice*
- Poster 36 Clare Pilgrim (MRC Clinical Sciences Centre, UK)
Nutrient sensing in the gut
- Poster 37 Joanna Poulton (University of Oxford, UK)
A common mtDNA variant predisposes to diabetes and impairs mitochondrial function. Studies in human mitochondrial cybrids, islets and muscle cultures
- Poster 38 Sara Ramírez (Universitat Internacional de Catalunya, Spain)
Hypothalamic ceramide levels regulated by CPT1C mediate orexigenic effect of ghrelin
- Poster 39 Cornelia Schöne (University of Cambridge, UK)
Fast and slow control of HA cells by optogenetic stimulation of hypocretin/orexin fibers
- Poster 40 Michelle Sim (University of Cambridge, UK)
Characterisation of pathogenic mutants of the lipodystrophy protein Seipin
- Poster 41 Hanna Sjoberg (University of Manchester, UK)
A systems-genetics approach to obesity and diabetes, using the BXD recombinant inbred model
- Poster 42 Lukasz Stasiak (University of Oxford, UK)
Functional analysis of fat mass and obesity associated (FTO) gene and protein
- Poster 43 Tatyana Tselovalnikova (Sechenov First Moscow State Medical University, Russia)
Metabolic abnormalities following treatment of brain tumors in young age
- Poster 44 Lourdes Valencia Torres (University of Cambridge, UK)
The role of 5-HT_{2C} receptors in the rewarding properties of food
- Poster 45 Amy Worth (University of Manchester, UK)
Postprandial effects of cholecystokinin are mediated by brainstem PrRP neurons
- Poster 46 Luigia Cristino (Italian Research Council, Italy)
Obesity-driven synaptic remodeling affects endocannabinoid control of orexinergic neurons