

Poster Index

- Poster 1 Christophe Blanchetot (arGEN-X BVBA, Belgium)
In vivo efficacy of human and mouse cross-reactive IL-22R SIMPLE antibody
- Poster 2 Emma Chambers (King's College London, UK)
IL-17A is increased in severe asthma and steroid refractory asthma: inhibition by 1alpha,25-dihydroxyvitamin D3 in CD39 dependent manner
- Poster 3 Susan Dallabrida (PHT Corporation, US)
Distribution of peak expiratory flow measurements from asthma patients captured wirelessly using an electronic handheld diary
- Poster 4 Nana Dang (Vrije Universiteit Brussel, Belgium)
Dendritic cells electroporated with IL-10 mRNA induce antigen-specific regulatory T cells
- Poster 5 Ismé de Kleer (Erasmus Medical Center, The Netherlands)
Highly increased IL33-ST2 axis in the lungs of neonatal mice increase susceptibility to house dust mite sensitization via CD11b+ dendritic cell skewing
- Poster 6 Lien Dejager (VIB, Ghent University, Belgium)
Neutralizing TNF restores glucocorticoid sensitivity in a mouse model of neutrophilic airway hyperinflammation
- Poster 7 Annelies Demeyer (VIB - DMBR, Belgium)
MALT1 deficiency in mice leads to regulatory T cell suppression with Th2 skewing, resulting in selective allergic skin pathology at later age
- Poster 8 Fien Devos (KU Leuven, Belgium)
Role of mast cells, IL-13 and TRP channels in a mouse model of chemical-induced airway hyperresponsiveness
- Poster 9 Anais Duval (CNRS, France)
IL-33 processing and activation by inflammatory proteases
- Poster 10 Gert Folkerts (University of Utrecht, The Netherlands)
Bifido Breve superior above Budesonide is a chronic model of allergic asthma?
- Poster 11 Carina Gross (NIBR Horsham, UK)
Impact of IL-13 on RSV infection in vitro and in vivo
- Poster 12 Rudi Hendriks (Erasmus MC Rotterdam, The Netherlands)
Enforced expression of Gata3 in the T cell and type 2 innate lymphocyte lineages increases susceptibility to allergic airway inflammation in mice
- Poster 13 Alex Kleinjan (Erasmus MC, The Netherlands)
Intervention of the notch pathway opens: new therapeutic entities in allergic asthma
- Poster 14 Mirjam Kool (Erasmus MC, The Netherlands)
Eosinophilic or neutrophilic asthma: a decisive role for the A20 enzyme in lung-specific dendritic cells
- Poster 15 Jisheng Liu (Ghent University, Belgium)
Permanent expression of BmToll9-1 receptor activates the innate immune response in silkworm-derived Bm5 cells

- Poster 16 Claire Mesnil (University of Leige, Belgium)
Resident CD11b+Ly6C-, not neo-recruited inflammatory CD11b+Ly6C+ dendritic cells mediate allergic airway sensitization to house dust mite antigens
- Poster 17 Mariel O'Connor (University of Nottingham, UK)
The effect of Azithromycin on proliferation and wound repair in human bronchial epithelial cells in vitro
- Poster 18 Zane Orinska (Research Center Borstel, Germany)
Modulation of IgE-mediated mast cell responses by anti-CD37 monoclonal antibodies
- Poster 19 Justyna Rzepecka (University of Strathclyde, UK)
From worm to pill: therapeutic effect of an ES-62-derived small molecule in experimental asthma
- Poster 20 Janeth Samson (Otsuka Pharmaceutical Inc., Philippines)
Effectiveness and safety of Budesonide Alone vs Budesonide/Formoterol in decreasing the number of severe exacerbations: a randomized controlled trial
- Poster 21 Marika Sarfati (CRCHUM, University of Montreal, Canada)
Basophils accumulated in inflamed mucosa tissues in Th2 as well as Th17/Th1 associated chronic inflammatory disorders in human and mice and directly promoted memory CD4 T cell responses
- Poster 22 Martijn Schuijs (VIB - DMBR, Belgium)
Role of GILZ in Th2 induction by DC's
- Poster 23 Sven Seys (KU Leuven, Belgium)
Sputum cytokine mapping reveals an "IL-5, IL-17A, IL-25-high" pattern associated with poorly controlled asthma
- Poster 24 Marina Shevchenko (Shemyakin and Ovchinnikov Institute of Bioorganic Chemistry, Russia)
The implication of extracellular Hsp70 in the natural regulation of allergic airway inflammation
- Poster 25 Anna Störmer (Research Center Borstel, Germany)
The allergy-protective Gram-positive cowshed bacteria Lactococcus lactis G121 and Staphylococcus sciuri W620 show differential mechanisms of immune cell activation and synergistic effects
- Poster 26 Cecilia, Mei Han Tong (University of Western Australia, Australia)
Immunomodulatory effects of the rye grass pollen Group 5 allergen on respiratory epithelial cells
- Poster 27 Heleen Vroman (Erasmus MC, The Netherlands)
Dendritic cell activation status directs T helper responses in a HDM allergic asthma model