

A model for accessibility and use of automated patch clamp system in academia: QPatch 48X in a core facility at UCL

University College London (UCL) - Sophion Consortium

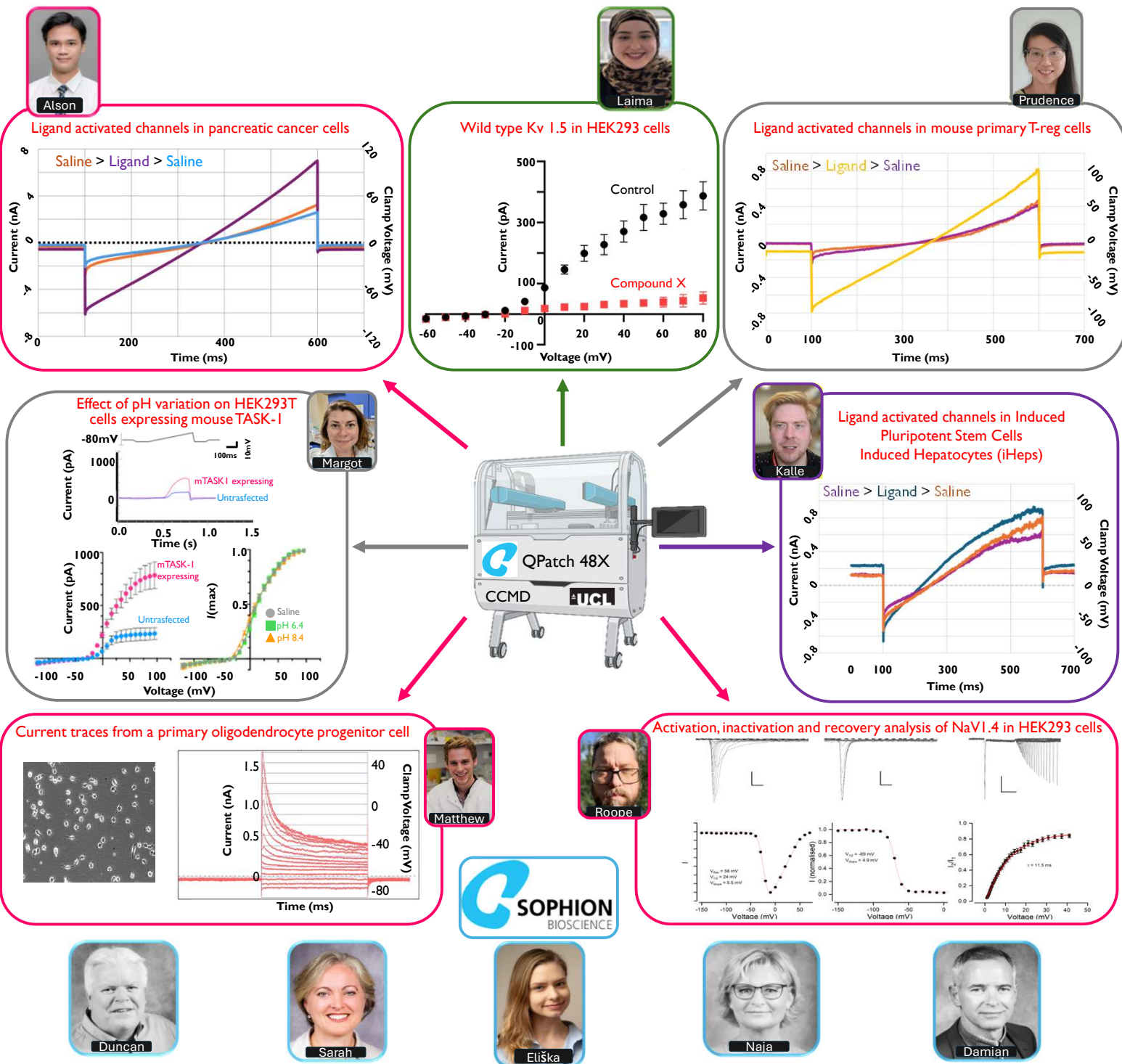
Centre for Cell and Molecular Dynamics, 5th Floor Rockefeller Building, University Street, University College London, London, UK

Sophion Biosciences A/S, Industriparken 39, 2750 Ballerup, Denmark



Aamir Ahmed*, Christopher Thrassivoulou*, Jonathan Ashmore*, Alson Huang*, Xiaohu Zhang*, Matthew Swire*, Arki Tian*, Stephanie Schorge*, Alasdair Gibb*, Dean Willis*, Ian Coombs*, Dorota Studniarczyk*, Roope Manniko*, Harry Luke-McClelland*, Conor McClenaghan*, Alan Greig*, Virginia Silo*, David Shorthouse*, Advay Ramgopal†, Kalle Sipilä†, Tamir Rashid†, Prudence Lui†, Niwa Ali†, Margot Maurer†, Mathilde Israel†, Stuart Bevan†, Kevin Cunningham†, Laima Tokhi†, Maryke Venter†, Duncan Jarman§, Sarah Lilley§, Naja Sørensen§, Damian Bell§, Eliška Waloschková§

To advance the use of automated patch clamp (APC) in academia, **Aamir Ahmed (UCL)** and **Sophion** joined forces to install a **QPatch 48X** in a core facility at the Centre for Cell and Molecular Dynamics (CCMD), UCL. The CCMD core facility allows easy access to QPatch for numerous local institutions with increased utility and less idle time for the machine. Since February 2025, with the support of colleagues from Sophion§, twelve different groups from four London universities (**UCL***, **Imperial College†**, **King's College‡** and **Westminster¶**) have used QPatch 48X for experiments with algae cells, primary mouse cells, human stem cells and cell lines. A selection of representative data are shown below.



We believe that a shared core facility model with multi-institutional user group, as established at UCL, provides a productive and cost-effective use of APC in academic institutions.