

# SEMINARS ON CELL MEMBRANE AND ION CHANNELS

## SEMINAR 1: SODIUM CHANNELS AND THE TUMOUR IONIC MICROENVIRONMENT

### INVITED SPEAKER

DR. WILLIAM J BRACKENBURY - UNIVERSITY OF YORK

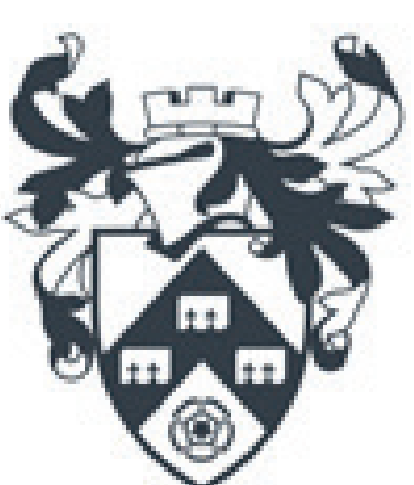
### MODERATOR

PROF. DR. GAMZE TANRIÖVER

DATE : 26 .09.2023  
TIME : 08.30-16.00  
VENUE : ISTANBUL MEDIPOL UNIVERSITY  
KAVACIK SOUTH CAMPUS  
CONFERENCE HALL



This event has been supported by TÜBİTAK (The Scientific and Technological Research Council of Turkey) within the scope of TÜBİTAK 2223-D Collaboration and Support for Organized Events Related to Priority Areas 2023/2 Call.



UNIVERSITY  
*of York*



MEDİPOL  
UNV-İSTANBUL  
İSTANBUL MEDİPOL ÜNİVERSİTESİ



# SEMINARS ON CELL MEMBRANE AND ION CHANNELS

## SEMINAR 2:

WELCOME TO THE BRAVE NEW\* WORLD OF  
AUTOMATED PATCH CLAMP:  
APC APPLICATIONS FURTHERING ION CHANNEL  
RESEARCH IN ACADEMIA & INDUSTRY

### INVITED SPEAKER

DR. DAMIAN BELL

### MODERATOR

DR. ŞERİFE YERLİKAYA

DATE : 26 .09.2023  
TIME : 08.30-16.00  
VENUE : ISTANBUL MEDIPOL UNIVERSITY  
KAVACIK SOUTH CAMPUS  
CONFERENCE HALL



This event has been supported by TÜBİTAK (The Scientific and Technological Research Council of Turkey) within the scope of TÜBİTAK 2223-D Collaboration and Support for Organized Events Related to Priority Areas 2023/2 Call.