

BPS19

63RD ANNUAL MEETING OF THE BIOPHYSICAL SOCIETY

BALTIMORE, MARYLAND • MARCH 2–6, 2019

nanjion

A NOVEL GAIN OF FUNCTION MUTATION OF PIEZO-1 IS INVESTIGATED IN RED BLOOD CELLS BY HIGH-THROUGHPUT PATCH CLAMP

Andrea Bruggemann, PhD¹, Giustina M. Rotordam¹, Nadine Becker, PhD¹, Niels Fertig, PhD¹, Paola Bianchi², Markus Rapedius, PhD¹, Lars Kaestner, PhD³.

¹Nanion Technologies, Munich, Germany, ²UOS Fisiopatologia delle Anemie, Fondazione IRCCS Ca' Granda Ospedale Maggiore Policlinico, UOC Ematologia, Milan, Italy, ³Theoretical Medicine and Biosciences, Homburg/Saar and Experimental Physics, Saarland University, Saarbrücken, Germany.

During their passage from the bone marrow to peripheral organs, Red Blood Cells (RBCs) experience substantial mechanical forces that trigger RBC volume changes via Ca²⁺ influx through mechanosensitive channels like Piezo-1 (1-3). However, direct electrophysiological proof of Piezo-1 activity in RBCs has been lacking so far as most of the studies come from animal models, were recorded under pathophysiological conditions or via Ca²⁺ imaging (1, 2, 4). In addition, the RBCs' small and variable size and their large heterogeneity in ion channel expression pose a significant challenge for successful patch clamp recordings (5, 7).

Here, we used automated patch clamp technology to study and characterize different ion channels present in RBCs, aiming specifically for Piezo-1 activity. We approached the problem of RBC heterogeneity by upscaling to high throughput automated patch clamp recordings of 384 cells in parallel to separate the cells in Yoda-1 responders and non-responders based on automated quality filtering. Thereby, we functionally characterized RBCs from patients carrying a gain of function mutation in Piezo-1 (8), demonstrating that high-throughput patch clamping can provide assays for drug discovery and personalized treatment of anaemic disorders such as hereditary xerocytosis and the sickle cell disease.

References:

- 1.) Danielczok et al., 2017
- 2.) Faucherre et al., 2014
- 3.) Gallagher et al., 2013
- 4.) Cahalan et al., 2015
- 5.) Bouyer et al., 2012
- 6.) Staines HM et al., 2007
- 7.) Minetti G et al., 2013
- 8.) Rotordam MG et al., 2018.