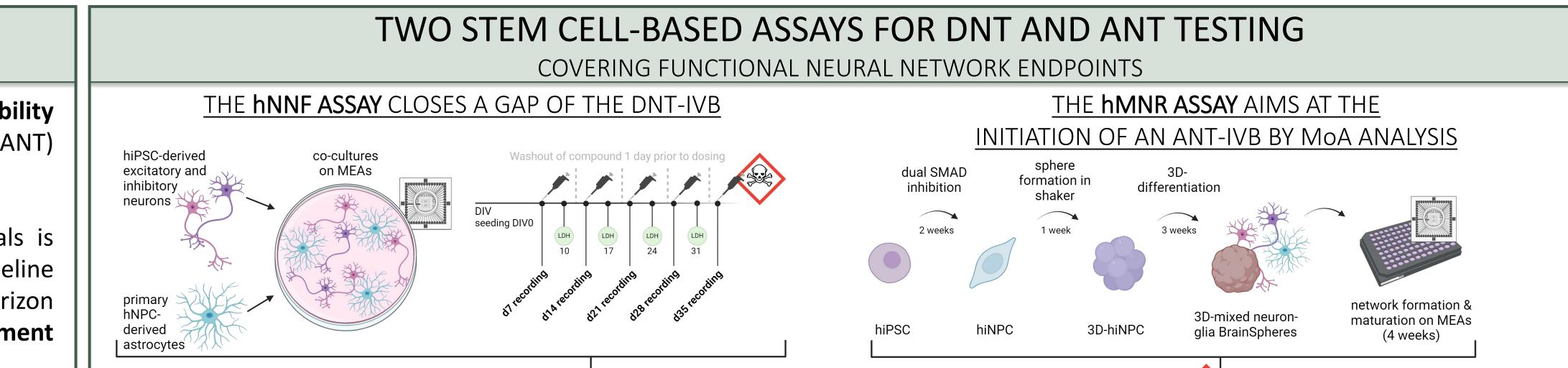


Refinement of stem cell-based in vitro assays towards a regulatory use for developmental and adult neurotoxicity testing of chemicals

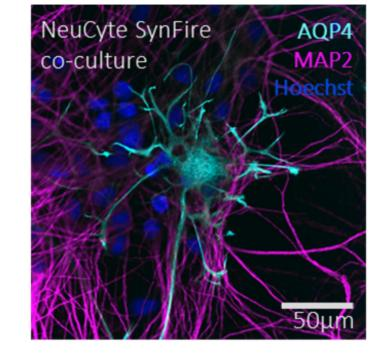
<u>Ilka Scharkin¹</u>, Kristina Bartmann^{1,2}, Sarah Delovec¹, Arif Dönmez^{1,2} Julia Hartmann¹, Etta Zühr¹, Volker Busskamp⁴, Andrea Rossi¹, *Katharina Koch^{1,2} and *Ellen Fritsche^{1,2,3} ¹ IUF - Leibniz Research Institute for Environmental Medicine, Duesseldorf, Germany; ³ Swiss Centre for Applied Human Toxicology (SCAHT), Basel, Switzerland; ⁴Department of Ophthalmology, Medical Faculty, University of Bonn, Bonn, Germany * Equal contribution

BACKGROUND & OBJECTIVES

- The EU Green Deal's Chemical Strategy for Sustainability highlights developmental (DNT) and adult neurotoxicity (ANT) as critical outcomes.
- Currently, regulatory DNT and ANT testing of chemicals is conducted in rats following OECD or US-EPA in vivo guideline studies. The **PARC Project** [1], initiated in 2022 under Horizon Europe, aims to implement Next Generation Risk Assessment using data from **New Approach Methods** (NAMs).



- An *in vitro* battery (DNT-IVB) has been established as an alternative for regulatory DNT testing [2]. However, neural network formation (NNF) in human-based systems was identified as a gap in the DNT-IVB v1.0, and there is currently no available IVB for ANT.
- To address this gap and to establish NAMs for an MoA-based first-generation ANT-IVB, two hiPSC-based assays have been developed: the human Neural Network Formation Assay (hNNF) for DNT and the human Multi-Neurotransmitter **Receptor Assay (hMNR)** for ANT, both using **microelectrode** arrays (MEAs).
- Both assays are undergoing **refinement** to reduce costs, increase throughput, and **improve regulatory acceptance**.



Assessment of spontaneous electrical activity Maestro Pro MEA System (Axion BioSystems)



Figure 1. The hNNF Assay. Commercially available, pre-differentiated hiPSC-derived excitatory and inhibitory neurons and primary human astrocytes (NeuCyte, USA) form functional co-cultures in 48-well (16 electrodes/well) MEA plates (Axion BioSystems). Cultures differentiate for 7 days in vitro (7DIV) before exposure to the test compound. 24h before the weekly recording of spontaneous network activity with the Maestro Pro MEA system (Axion BioSystems), a washout of the compound is performed. Together with a weekly cytotoxicity assessment, this procedure is repeated until DIV35. Corresponding publication: [3] Bartmann et al. (2023).

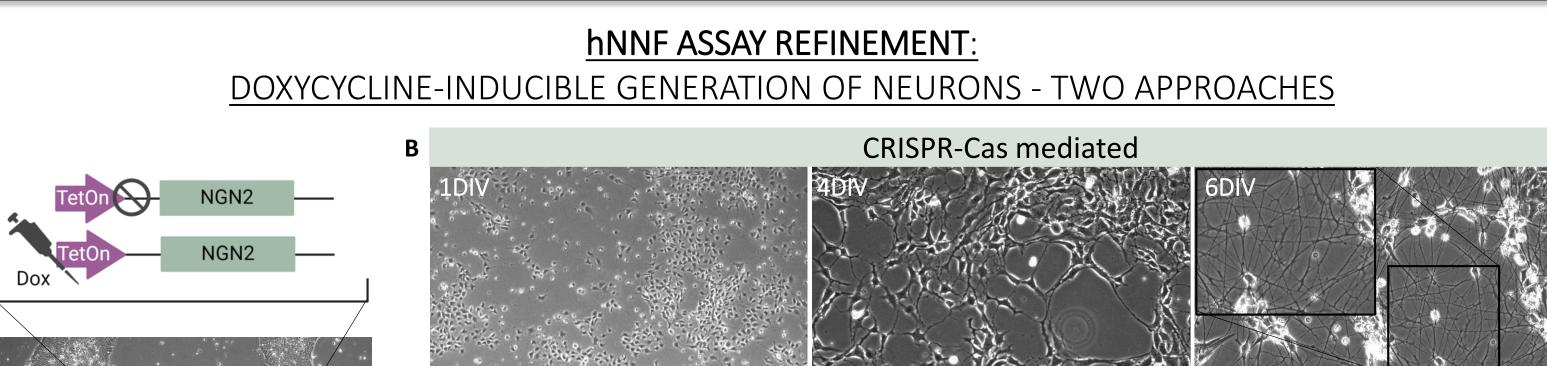
Assessment of spontaneou electrical activity Maestro Pro MEA System (AxionBioSystems)

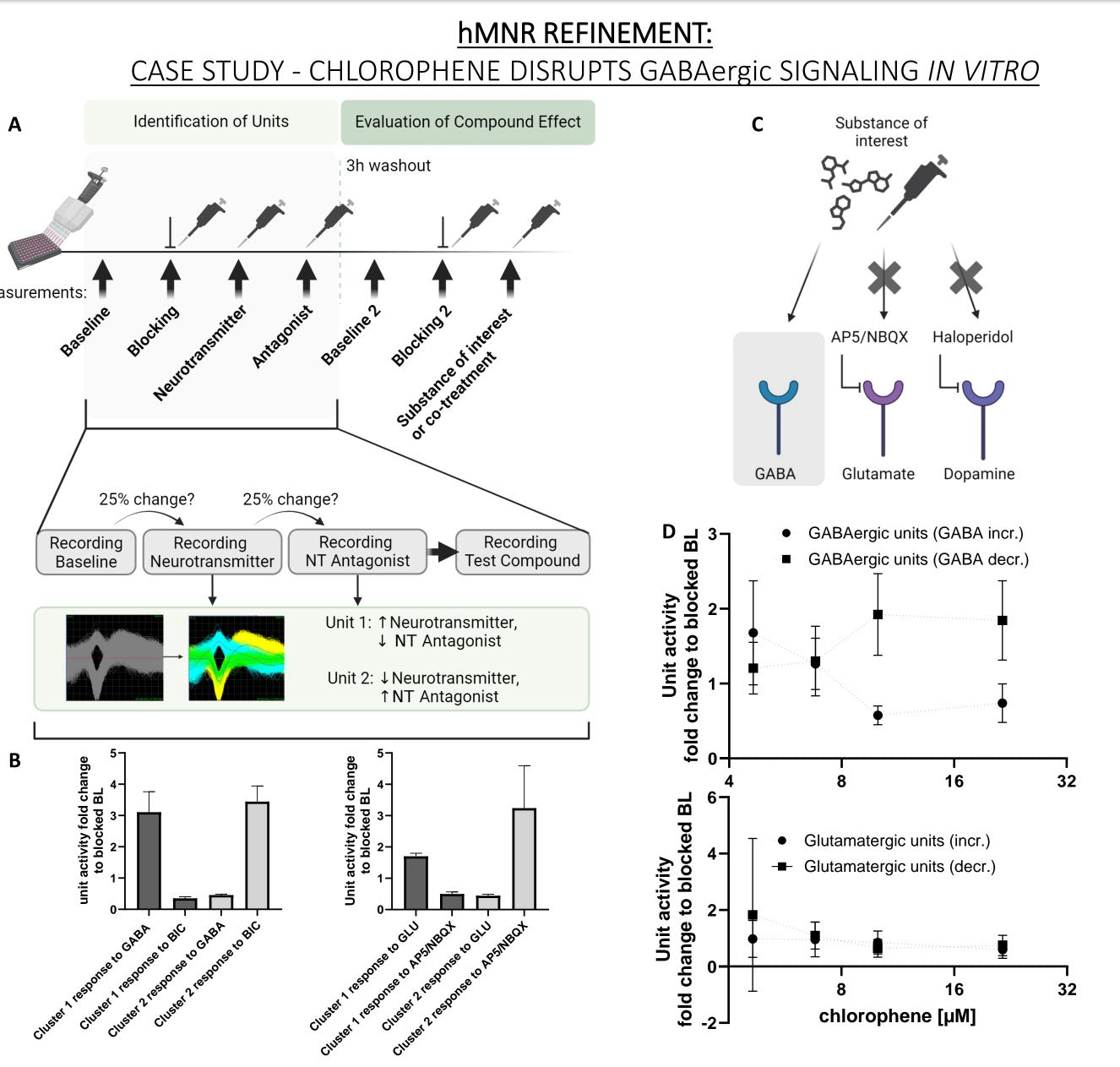
Figure 2. The hMNR Assay. Dual SMAD inhibition of hiPSCs generates hiNPCs. Shaking cultivation in CINDA+ differentiation medium yields 3D mixed neuron-glia BrainSpheres containing glutamatergic, GABA-ergic, dopaminergic, serotonergic and cholinergic neuronal subtypes that exhibit spontaneous electrical activity on 96-well MEAs. Detected spikes are sorted using the Plexon Offline Sorter software after treatment with neurotransmitter receptor agonists/antagonists. Compound effects on neuronal subtype receptors, including GABAergic, glutamatergic, and dopaminergic units, can be detected and quantified. Corresponding publication: [4] Hartmann et al. (2023).

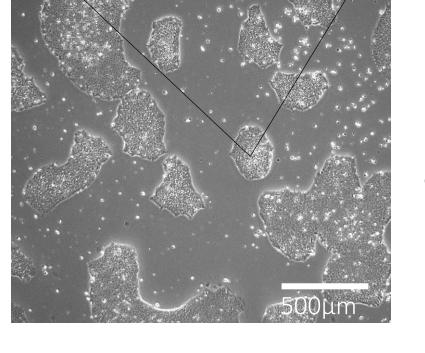
Abbreviations: hiPSC - human induced pluripotent stem cell; hNPC - human neural progenitor cells; MEA - microelectrode array; Prol. - proliferation; S100ß - S100 calcium binding protein ß (Astrocyte marker); TUBB3 - *β*-III-tubulin (Neuron marker); AQP4 - Aquaporin 4 (Astrocyte marker), MAP2 - Microtubule-associated protein 2 (Neuron marker)

RESULTS

ASSAY REFINEMENT TOWARDS ENHANCED THROUGHPUT AND COST REDUCTION







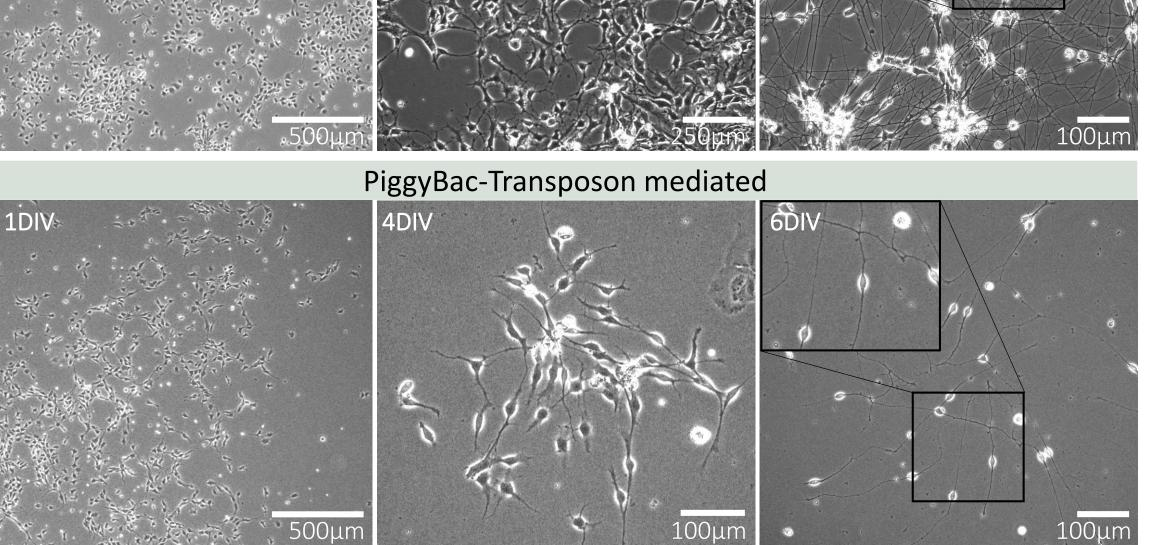


Figure 3. In-house-generated hiPSC lines differentiate into post-mitotic excitatory (Ngn1/2) neurons through doxycycline induction. Two approaches for generating inducible neurons are being tested: Commercially available IMR90 hiPSCs (A, WiCell, USA) are equipped with a doxycycline-inducible expression cassette for the excitatory neuronal transcription factor Ngn1/2 by CRISPR-Cas (B) or the PiggyBac-Transposon System (C). With both methods, post-mitotic, network-forming neuronal cultures have been generated within six days in our laboratories. *Corresponding publication:* [5] *Busskamp et al. (2014).*

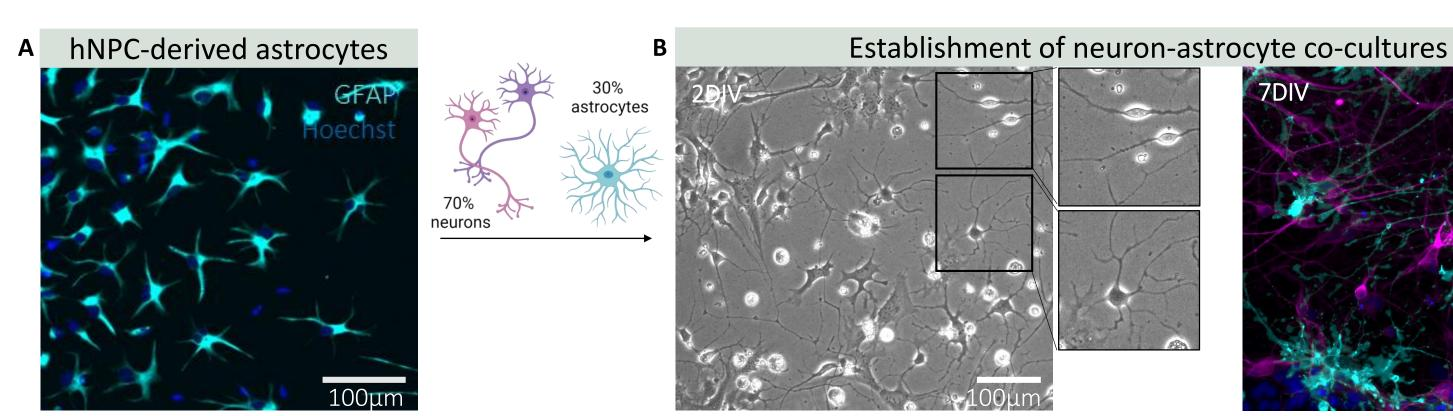


Figure 4. Neuron-astrocyte co-cultures. Combined with a fast hNPC-based astrocyte differentiation protocol (A, established by Etta Zühr, IUF), first successful defined co-cultures using the CRISPR-Cas-generated excitatory neurons (see Fig. 3B) have been established. Pre-differentiated hiPSC-based neurons together with cryopreserved and batched hNPC-based astrocytes were co-cultured in a 70%/30% ratio. (B) After two days of culture, both cell types present the characteristic morphologies (see enlarged boxes: top – neurons, bottom: astrocytes). Performing immunocytochemistry at day 7 in culture, both neurons and astrocytes could be identified by the corresponding markers MAP2 and GFAP.

Figure 5. hMNR case study: Chlorophene disrupts GABAergic signaling in vitro using the refined hMNR-Assay. (A) The assay workflow involves spike sorting with the Plexon Offline Sorter (Plexon, USA) software to identify units reacting to receptor agonists and antagonists prior to compound effect analysis. A cutoff of 25% change in spiking between baseline and neurotransmitter addition determines unit identification. (B) This approach can discriminate early and later stages of neuronal subtype maturation, like pre- and post-GABA-switch neurons. Displayed as mean +- SEM. n >= 21 (units) per response cluster. (C) Refinements include an additional blocking step for precise measurement of signals from the neuronal subtype of interest, exemplified for GABAergic units. (D) Collaboration with the UFZ Leipzig using BrainSpheres treated with Chlorophene in the hMNR Assay (manuscript in prep.) reveals opposing responses in pre- and post-GABA switch-like units, with glutamatergic units unaffected. Displayed as mean +- SEM . $n \ge 4$ (units) per response cluster. Ongoing refinement aims to develop a robust data evaluation strategy for higher-throughput analysis of MEA data.

Abbreviations: DIV - days in vitro; Ngn1/2 - Neurogenin1/2; CRISPR-Cas - Clustered regulalrly interspaced short palindromic repeats and CRISPR-associated protein; GFAP - glial fibrillary acidic protein; MAP2 - microtubule-associated protein; BIC - Bicuculline, GABA A receptor antagonist; AP5/NBQX - glutamate receptor antagonists; GLU - glutamate, incr. - increase; decr. - decrease.

CONCLUSION & OUTLOOK	LINKS & LITERATURE
 Set-up of the hNNF assay closed one identified gap of the DNT-IVB. The ongoing refinement using in-house generated hiPSC-based cell lines aims at promote cost-efficient alternative, enabling a higher throughput and extensive test system characterization. 	roviding a [1] Marx-Stoelting, P. et al. (2023) doi: 10.1007/s00204-022-03435-7. [2] Blum, J. et al. (2022)
• There is a need for setting up a first-generation ANT-IVB with test methods focusing on known MoA of neurotoxicity. The refined hMNR assay aims efficient and unified evaluation of large data sets from functional MEA readouts, inclusion of male and female donor-derived hiPSC lines as well a characterization of the cell system for more exact determination of applicability domains. Testing of model compounds demonstrated the suitability of the the MoA identification for multiple neurotransmitter receptors [3].	as further doi: 10.14573/altex.2206031. (4] Hartmann, J. et al. (2023) doi: 10.3390/cells12091270. (5] Busskamp, V. et al. (2014)
• The described in vitro test methods will further contribute to the establishment and acceptance of DNT- and ANT-IVBs. Efforts continue to refine bot towards regulatory suitability and acceptance, with future validation planned through testing of known positive and negative DNT and ANT model compositive and negative DNT and	
towards regulatory suitability and acceptance, with future validation planned through testing of known positive and negative DNT and ANT model compounds. All figures created with biorender.com. Contact: Ilka Scharkin (Ilka.Scharkin@iuf-duesseldorf.de) IUF – Leibniz Research Institute for Environmental Medicine Image: Contact in the second	

This work was funded by Horizon Europe PARC (Grant Agreement No. 101057014), the Danish EPA (Grant No. MST-66-00205) and CERST Germany (File No. 233-1.08.03.03-121972).