

Effects of mexiletine and its pyrroline derivative on Na_v1.4 and Na_v1.5 by automated patch-clamp: towards safer anti-myotonic drugs

Alessandro Giovanni Cerchiara¹, Nadine Becker², Niels Fertig², Ornella Cappellari¹, George Okeyo³, Michela De Bellis¹, Alessia Carocci¹, Giovanni Lentini¹, Jean-Francois Rolland⁴, Paola Imbrici¹, Annamaria De Luca¹, Alison Obergrussberger²

¹ Department of Pharmacy – Drug Sciences, University of Bari "Aldo Moro", Bari, Italy; ² Nanion Technologies GmbH, Munich, Germany; ³ Nanion Technologies Inc. Livingston, NJ, USA; ⁴ Axxam, SpA, Bresso Italy

Contact: ali@nanion.de, annamaria.deluca@uniba.it;

nanjion

UNIVERSITÀ
DEGLI STUDI DI BARI
ALDO MORO



1. Introduction

Myotonia is an inherited rare skeletal muscle disorder which causes sarcolemma hyperexcitability due to alterations of different ion channels (e.g., CLC-1, Na_v1.4)¹. Mexiletine is a class I antiarrhythmic drug approved in non-dystrophic myotonias for its ability to exert a use-dependent block of skeletal muscle sodium channel (Na_v1.4)^{2,3}. Mexiletine showed to reduce muscle stiffness, weakness and pain in patients of two randomized controlled clinical trials^{4,5} and to be safe in two long-term clinical studies^{6,7}. Nevertheless, due to different side effects (e.g. gastrointestinal and cardiac problems, headache, dizziness, etc.) or lack of response, several patients are not eligible for mexiletine treatment^{4,6}. Consequently, drug potency towards Na_v isoforms remains an open issue for tissue selectivity and efficacy/safety ratio. Moreover, preclinical studies showed that other antiarrhythmic drugs or their derivatives presented improved efficacy in myotonic conditions, compared to mexiletine^{8,9,10}.

The possibility to discriminate drug action between Na_v1.4 and cardiac Na_v1.5 is critical for exerting a therapeutic anti-myotonic effect, reducing the risk of cardiac toxicity. Efficient and reliable pre-clinical tests are pivotal, also with the long-term goal to find novel anti-myotonic agents. Furthermore, use of alternative drugs to safely treat myotonia could be pivotal for more efficient therapy. Our aim was to assess effects mexiletine and its derivative compound (VM11) both on Na_v1.4 and Na_v1.5 using an automated patch clamp platform.

2. Methods

Cell culture

We used TE671 cells (provided by Cell Line Services: <https://cls.shop/>), which endogenously express Na_v1.4, and CHO cells heterologously expressing Na_v1.5 (provided by Charles River: <https://www.criver.com/find-cell-line>).

Compounds

The compounds tested were 2-(2,6-dimethylphenoxy)-1-amino-propanemethylethylamine (Mex) and the mexiletine-derivative 2,2,5,5-tetramethyl-N-[1-methyl-2-(2,6-dimethylphenoxy)ethyl]-2,5-dihydro-1H (VM11). Stock solutions containing external solution and dimethyl sulfoxide (<0.33%) were used for both compounds.

Electrophysiology

Recordings of peak sodium currents (at -20 mV) of Na_v1.4 and Na_v1.5 were performed using the automated patch clamp platform Patchliner. For dose response curves, at least 8 increasing concentrations were included. The tonic (0.3 Hz), voltage- (-70 and -140 mV at 0.3 Hz) and use-dependent blocks (3 Hz for Na_v1.5 and 10 Hz for Na_v1.4), consistent with pathophysiological conditions of Na_v1.4 and Na_v1.5 by Mex/VM11, were evaluated.



Figure 1. Patchliner: fully automated planar patch clamp system

3. Tonic (TB) and voltage-dependent (VDB) blocks

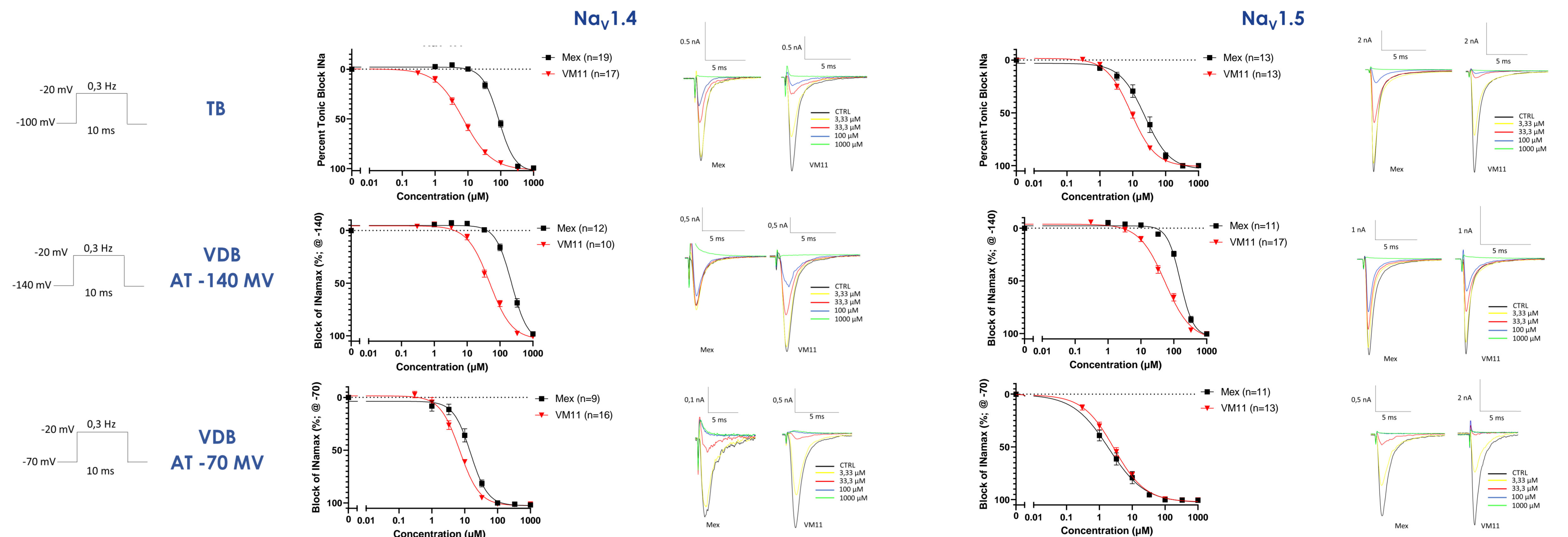


Figure 2: Dose-response curves depicting the tonic (A) and voltage-dependent block at -70 (B) and -140 mV (C) of Mex and VM11 for Na_v1.4 and 1.5. Mex was more potent in determining a TB of Na_v1.5 than of Na_v1.4 measured at 0.3 Hz (IC₅₀: 7 ± 0.6 vs 87 ± 5 μM). VDB at -140 mV showed that Mex was more potent in blocking Na_v1.5 than Na_v1.4. (159 ± 8 vs 226 ± 16 μM). On the other hand, VM11 produced similar TB and VDB (-140mV) of Na_v1.4 and Na_v1.5 (7 ± 0.6 vs 9 ± 0.5 μM for TB; 48 ± 3 vs 54 ± 5 μM for VDB). Mex produced a more potent VDB (at -70mV) on Na_v1.5 than Na_v1.4 (2 ± 0.3 vs 15 ± 1 μM). Similarly, VM11 blocked Na_v1.5 more than Na_v1.4 in VDB (at -70mV) conditions (3 ± 0.2 vs 7 ± 0.4 μM).

4. Use-dependent block (UDB)

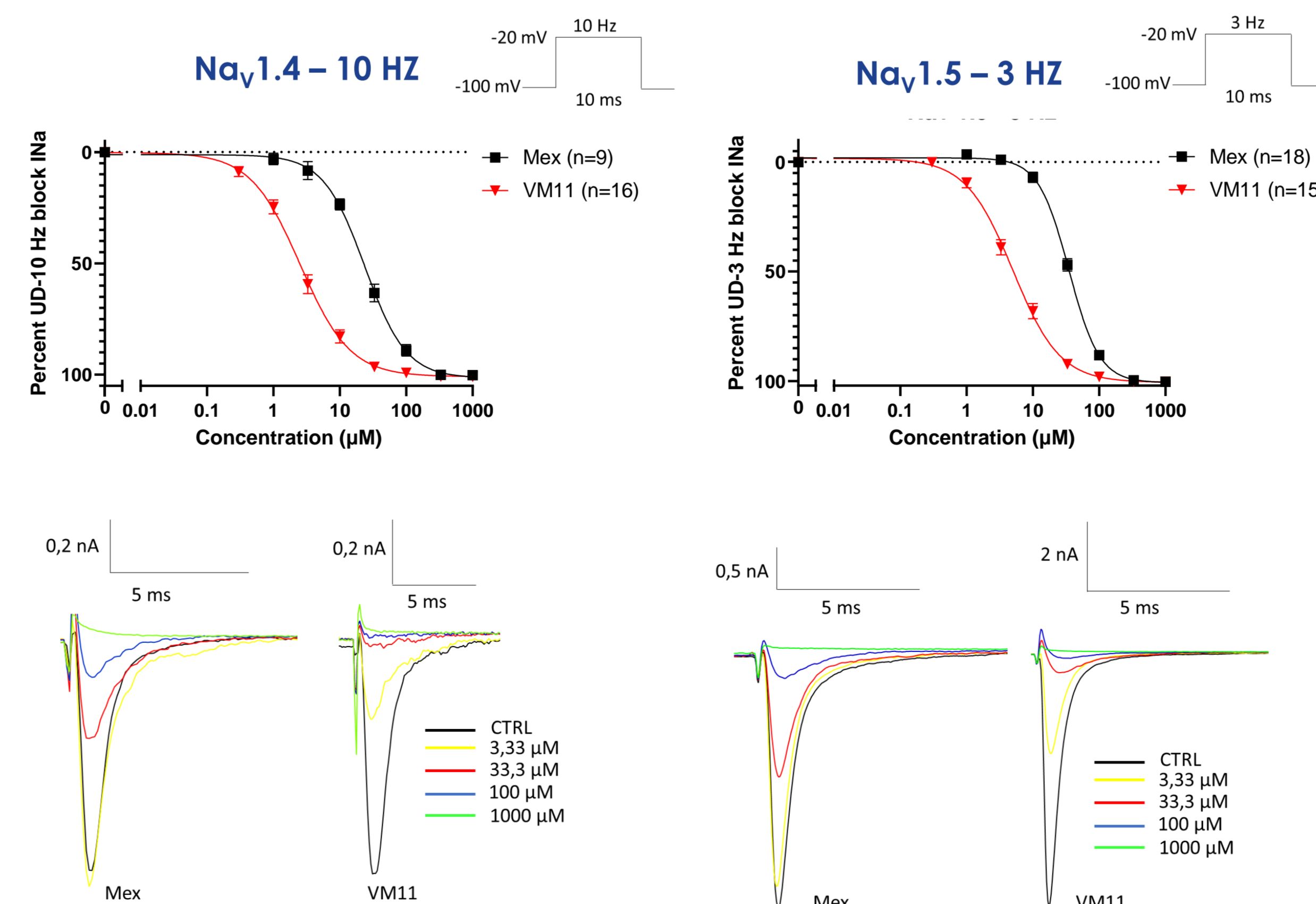


Figure 3. Dose-Response curves showing UDB of Mex and VM11 on Na_v1.4 and Na_v1.5.

At cardiac pathophysiological frequency of 3 Hz, VM11 blocked Na_v1.5 more than Mex with an IC₅₀ of 5 ± 0.3 μM (vs 35 ± 1 μM). At stimulation frequencies mimicking the myotonic discharge (10 Hz), both Mex and VM11 blocked Na_v1.4 channels with high potency (24 ± 2 and 3 ± 0.2 μM).

5. Conclusions

- ✓ Mex has more affinity for Na_v1.5 channel compared to Na_v1.4 in conditions of TB/VDB.
- ✓ VM11 was more potent in determining TB and VDB (at -140 mV) of Na_v1.4 than Na_v1.5, but blocks Na_v1.5 more than Na_v1.4 in condition of VDB at -70 mV.
- ✓ At pathological heart frequencies (3 Hz), VM11 blocked Na_v1.5 more than Mex, while at myotonic-like frequency of 10 Hz, Mex and VM11 potently blocked Na_v1.4 but still at concentrations in the range of blocking Na_v1.5.
- ✓ This biophysical approach can help the development of novel Na_v1.4 blockers with improved safety profile for muscle disorders.

References

1. Stunnenberg et al. *Muscle & nerve*. 2020;62(4):430-444.
2. De Luca et al. *Neuromuscular disorders*. 2004;14(7):405-416.
3. Hoffman & Kaminski. *JAMA*. 2012;308(13):1377-1378.
4. Statland et al. *JAMA*. 2012;308(13):1357-1365.
5. Stunnenberg et al. *JAMA*. 2018;320(22):2344-2353.
6. Suetterlin et al. *JAMA neurology*. 2015;72(12):1531-1533.
7. Modoni et al. *Frontiers in neurology*. 2020;11:300.
8. Desaphy et al. *The Journal of physiology*. 2004;554(Pt 2):321-334.
9. De Bellis et al. *Neuropharmacology*. 2017;113(Pt A):206-216.
10. De Bellis et al. *Frontiers in pharmacology*. 2017;8:907.

