



Buffer solutions

Designed to set new standards



nanjion

Key benefits at a glance

- | | | | |
|----------|--|----------|--|
| 1 | Tested high quality solutions with a long shelf life | 5 | ISO-certified, GMP-compliant, certified analysis and detailed MSDS |
| 2 | No batch-to-batch variability | 6 | Highest chemical purity on the market (ACS, ISO, Ph. Eur.) |
| 3 | High throughput APC without fluoride | 7 | Tailored bulk purchase discounts available |
| 4 | Ready-to-use, convenient packaging, storage at RT | 8 | Recyclable PET-bottles |



High quality reagents for automated patch clamp

Quality-assured and ISO-certified

Innovations in electrophysiology have prompted a high demand for simplified workflows and solutions you can rely on. Nanion's stable, long-life, and ready-to-use buffer solutions are produced to the highest quality under ISO-certified and GMP-compliant conditions, ensuring reliability and excellence in your research.

Increase your productivity

Nanion's buffer solutions are certified, tested and made with the highest purity chemicals. Our wide range of solutions, short delivery times and tailored discount schemes ensure that you meet your research goals ontime and within budget. All our solutions are quality-checked at our headquarters and provided in recyclable PET bottles.

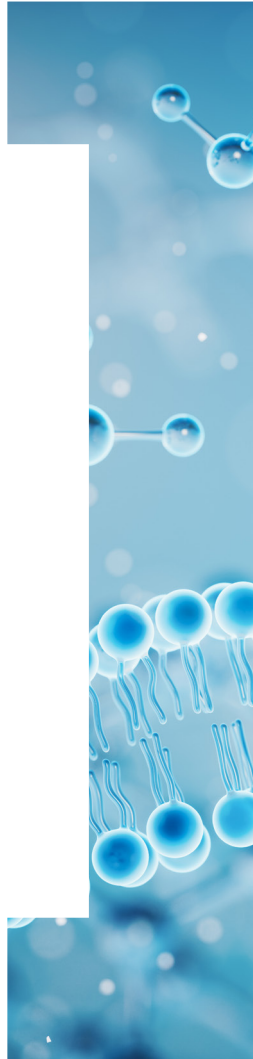


Our solutions are manufactured in cooperation with award-winning and certified production facilities (PAN-Biotech GmbH, Germany and Ampliqon A/S, Denmark).

Solution	Color code	Order #	Description/Based on
External Standard	●	08 3001	NaCl, Hepes
External Standard Ca 10	●	08 3012	NaCl, Ca ²⁺ , Hepes
External [-] Ca ²⁺ [-] Mg ²⁺	●	08 3003	Divalent free
External NMDG 60	●	08 3004	NMDG
External NMDG 60 Ca 10	●	08 3011	NMDG, Ca ²⁺
External NMDG 140	●	08 3014	For flouride free measurements
Washing Solution for Tips	●	08 3010	24 % EtOH, EDTA
Internal CsF 110	●	08 3008	High CsF, EGTA
Internal KF 110	●	08 3007	High KF, EGTA
Internal K Gluconate 120	●	08 3015	For flouride free measurements
Reagent Kit Port-a-Patch standard	●	08 1001	Starter Kit with 81 vials
Reagent Kit Port-a-Patch for Na _v	●	08 1014	Starter Kit with 81 vials



accelerate your research



nanion Europe

info@nanion.de

phone: +49 89 2190 95-0

www.nanion.de

nanion USA

info@naniontech.com

phone: 1-888-9-NANION

www.nanion.de

nanion China

andy.di@nanion.cn

phone: +86 10 82 17 6388

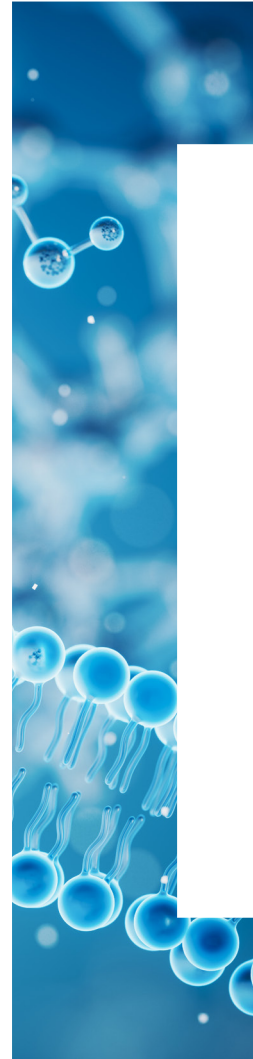
www.nanion.de/zh-hans

nanion Japan

info@nanion.jp

phone: +81 3 6555 3773

www.nanion.de/ja



To find your nearest Nanion
representative, visit our website:
www.nanion.de

nan]i[on