

Product Specification:



QCells

CHO-Na_v1.5 optimized for the QPatch



- Cells optimized for the QPatch
- Tight sealing properties
- Optimal success rate
- Stable current response
- Cell line support
- Optimized Ringer's solutions
- Custom assay and analysis

See specifications on back

The sodium ion channel $\text{Na}_v1.5$ is expressed as an integral membrane protein and contains a tetrodotoxin-resistant voltage-gated sodium channel subunit.

The encoded protein is found primarily in cardiac muscle and is responsible for the initial upstroke of the action potential in an electrocardiogram

Mutations in the gene are associated with long QT syndrome type 3, Brugada syndrome, primary cardiac conduction disease and idiopathic ventricular fibrillation.

