

# Product Information



## Ca<sub>v</sub>1.2 Calcium Channel Monoclonal Antibody (Clone S57-46)

Catalog No. 13702

<b>Contents:</b>	This vial contains 100 µg of protein G-purified IgG in 100 µl PBS, pH 7.4, containing 50% glycerol and 0.09% sodium azide.
<b>Antigen:</b>	Fusion protein from rabbit Ca <sub>v</sub> 1.2 amino acids 1,507-1,733
<b>Isotype:</b>	IgG <sub>2b</sub>
<b>Host:</b>	Mouse, clone S57-46
<b>Cross Reactivity:</b>	(+) Human, murine, and rat Ca <sub>v</sub> 1.2
<b>Stability:</b>	≥1 year at -20°C
<b>Applications:</b>	Western blot (WB), immunohistochemistry (IHC), and immunocytochemistry (ICC). The recommended starting dilution for WB is 1-10 µg/ml, IHC/ICC is 0.1-1.0 µg/ml (HRP detection), and IF is 1-10 µg/ml.

Ion channels are integral membrane proteins that help establish and control the small voltage gradient across the plasma membrane of living cells by allowing the flow of ions down their electrochemical gradient.<sup>1</sup> They are present in the membranes that surround all biological cells because their main function is to regulate the flow of ions across this membrane. Whereas some ion channels permit the passage of ions based on charge, others conduct based on an ionic species, such as sodium or potassium. Furthermore, in some ion channels, the passage is governed by a gate which is controlled by chemical or electrical signals, temperature, or mechanical forces.

There are a few main classifications of gated ion channels. There are voltage-gated ion channels, ligand-gated, other gating systems, and finally those that are classified differently, having more exotic characteristics. The first are voltage-gated ion channels which open and close in response to membrane potential. These are then separated into sodium, calcium, potassium, proton, transient receptor, and cyclic nucleotide-gated channels, each of which is responsible for a unique role. Ligand-gated ion channels are also known as ionotropic receptors, and they open in response to specific ligand molecules binding to the extracellular domain of the receptor protein. The other gated classifications include activation and inactivation by second messengers, inward-rectifier potassium channels, calcium-activated potassium channels, two-pore-domain potassium channels, light-gated channels, mechano-sensitive ion channels, and cyclic nucleotide-gated channels. Finally, the other classifications are based on less normal characteristics such as two-pore channels, and transient receptor potential channels.<sup>2</sup>

Specifically, Ca<sub>v</sub>1.2 is a cardiac L-type calcium channel and is important for excitation and contraction of the heart.<sup>3</sup> It may be associated with a variant of Long QT syndrome called Timothy's syndrome<sup>4,5</sup> and also with Brugada syndrome. Some references suggest it is related to bipolar disease as well.<sup>5</sup>

### References

- Hille, B. Ion Channels of Excitable Membranes. 3rd Ed., Sinauer Associates Inc., Sunderland, MA (2001).
- What are ion channels? Retrieved October 22, 2009, from <http://www.ionchannels.org/>.
- Splawski, I., Timothy, K.W., Sharpe, L.M., *et al.* Ca<sub>v</sub>1.2 calcium channel dysfunction causes a multisystem disorder including arrhythmia and autism. *Cell* **119**, 19-31 (2004).
- Krey, J.F., Dolmetsch, R. The Timothy Syndrome mutation in Ca<sub>v</sub>1.2 causes dendritic retraction through calcium-independent activation of the rho A pathway. *Biophysical* **96(3)**, 221a-222a (2009).
- Crotti, L., Celano, G., Dagradi, F., *et al.* Congenital long QT syndrome. *Orphanet Journal of Rare Diseases* **3**, 18-31 (2008).

### Related Products

Ca<sub>v</sub>β1 Calcium Channel Monoclonal Antibody (Clone S7-18) - Cat. No. 13700 • Ca<sub>v</sub>β4 Calcium Channel Monoclonal Antibody (Clone S10-7) - Cat. No. 13701 • Ca<sub>v</sub>1.3 Calcium Channel Monoclonal Antibody (Clone S38-8) - Cat. No. 13703 • Ca<sub>v</sub>3.2 Calcium Channel Monoclonal Antibody (Clone S55-10) - Cat. No. 13704 • Ca<sub>v</sub>1.3 Calcium Channel Monoclonal Antibody (Clone S48A-9) - Cat. No. 13706

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 11/20/2009

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)