

PRODUCT INFORMATION

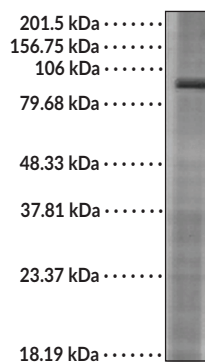
HCN2 Cyclic Nucleotide-gated Channel Monoclonal Antibody (Clone S71-37)

Item No. 13707

Overview and Properties

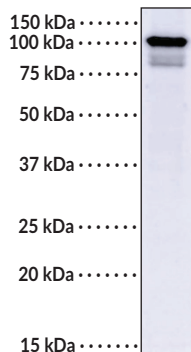
Contents:	This vial contains 100 µg of protein G-purified monoclonal antibody.
Immunogen:	Fusion protein amino acids 761-863 (cytoplasmic C-terminus) of rat HCN2
Species Reactivity:	(+) Human, mouse, rat; other species not tested
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	100 µl PBS, pH 7.4, containing 50% glycerol and 0.09% sodium azide
Concentration:	1 mg/ml
Clone:	S70
Host:	Mouse
Isotype:	IgG1
Applications:	Immunocytochemistry (ICC), Immunohistochemistry (IHC), Immunoprecipitation (IP), and Western blot (WB); the recommended starting dilution is 1:100 for ICC and 1:1,000 for IHC and WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: HCN2 Cyclic Nucleotide-gated Channel (15 µg)

WB of HCN2 Cyclic Nucleotide-gated Channel Monoclonal Antibody (Clone S71-37).



Lane 1: HCN2 Cyclic Nucleotide-gated Channel (15 µg)

WB of mouse brain showing detection of approximately 95 kDa HCN2 protein using HCN2 Cyclic Nucleotide-gated Channel Monoclonal Antibody (Clone S71-37).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/30/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

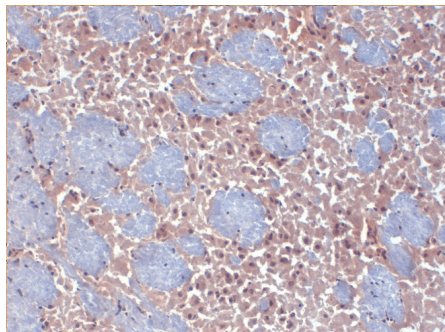
CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

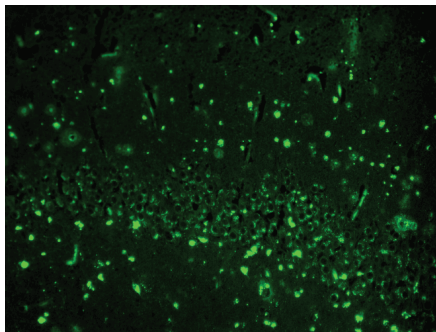
PRODUCT INFORMATION



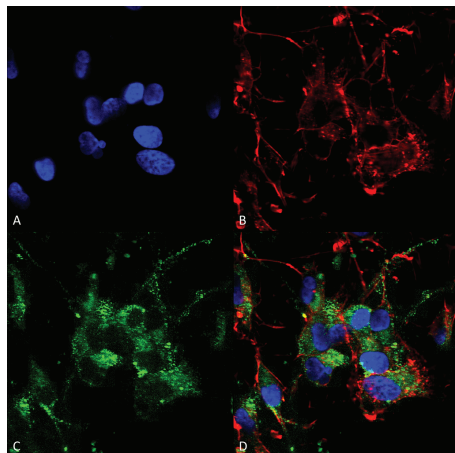
Images



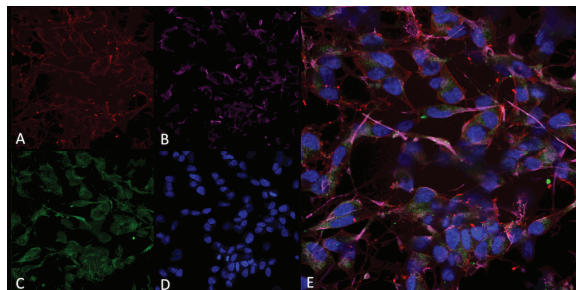
Immunohistochemical staining of formalin-fixed mouse brain using HCN2 Cyclic Nucleotide-gated Channel Monoclonal Antibody (Clone S71-37) at a dilution of 1:1,000 for one hour at room temperature. Following incubation with HCN2 Cyclic Nucleotide-gated Channel Monoclonal Antibody (Clone S71-37), cells/tissues were incubated with biotinylated goat anti-mouse, streptavidin peroxidase, DAB chromogen (brown) for 30 minutes at room temperature and stained with Mayer Hematoxylin (purple/blue) nuclear stain at 250-500 μ l for five minutes at room temperature.



Immunohistochemical staining of Bouin's fixative and paraffin-embedded human hippocampus using HCN2 Cyclic Nucleotide-gated Channel Monoclonal Antibody (Clone S71-37) at a dilution of 1:100 for one hour at room temperature. Following incubation with HCN2 Cyclic Nucleotide-gated Channel Monoclonal Antibody (Clone S71-37), tissues were incubated with FITC goat anti-mouse (green) for one hour at room temperature.



Immunocytochemical/Immunofluorescence staining of formalin-fixed human neuroblast using HCN2 Cyclic Nucleotide-gated Channel Monoclonal Antibody (Clone S71-37) at a dilution of 1:50 for overnight at 4°C with slow rocking. Following incubation with HCN2 Cyclic Nucleotide-gated Channel Monoclonal Antibody (Clone S71-37), cells/tissues were incubated with AlexaFluor 488 at a dilution of 1:1,000 for 1 hour at room temperature and stained with Phalloidin-iFluor 647 (red) F-Actin stain, Hoechst (blue) nuclear stain at 1:800, 1.6 mM for 20 minutes at room temperature with the following: (A) Hoechst (blue) nuclear stain (B) Phalloidin-iFluor 647 (red) F-Actin stain (C) HCN1 Antibody (D) Composite.



Immunocytochemical/Immunofluorescence staining of formalin-fixed human Differentiated SH-SY5Y using HCN2 Cyclic Nucleotide-gated Channel Monoclonal Antibody (Clone S71-37) at a dilution of 1:100. Following incubation with HCN2 Cyclic Nucleotide-gated Channel Monoclonal Antibody (Clone S71-37), cells/tissues were incubated with AlexaFluor 488 and stained with Phalloidin-iFluor 647 (red) F-Actin stain, Hoechst (blue) nuclear stain with the following: (A) Phalloidin (B) Anti-beta III Tubulin Ab. (C) HCN2 Antibody (D) Hoechst (E) Composite.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/30/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Hyperpolarization-activated cation channels of the HCN gene family contribute to spontaneous rhythmic activity in both the heart and the brain.¹

Reference

1. Zong, X., Eckert, C., Yuan, H., *et al.* A novel mechanism of modulation of hyperpolarization-activated cyclic nucleotide-gated channels by Src kinase. *J. Biol. Chem.* **280**(40), 34224-34232 (2005).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM