

# PRODUCT DATA SHEET



**CATALOG NO.:** SA-598

**LOT NO.:** temp

**PRODUCT:** Anti-HCN2 (Hyperpolarization-activated Cyclic Nucleotide-Gated Potassium Channel 2, HAC1, BCNG2) and Control Peptide.

KI-425 Anti-HCN2

KI-426 HCN2 Control Peptide, 40 µg

**IMMUNOGEN:** A synthetic peptide corresponding to amino acid residues 147-161 of human HCN2, Accession: CAB42602. This sequence is identical in mouse and rat HCN2.

**SPECIFICITY:** Recognizes HCN2 from rat and human tissue.

**APPLICATIONS:** WB: 1:200-1000, IH.

**FORM / PURIFICATION:** Rabbit polyclonal antibody: Affinity purified IgG. Supplied lyophilized from PBS, Ph 7.4, 1% BSA and 0.05% sodium azide. After reconstitution antibody concentration is **0.8 mg/ml**. Control peptide: Supplied lyophilized.

**RECONSTITUTION:** To lyophilized antibody, add 50 µl or 200 µl (depending on which size was purchased) of deionized water. To lyophilized control peptide, add 100 µl of distilled water.

**PREADSORPTION CONTROL:** 1 µg control peptide per 1 µg antibody. Prepare 2 tubes. In the first, dilute the necessary amount of antibody in 200-500 µl of PBS, 1% BSA and 0.025% sodium azide. Prepare the second tube identically to the first with the addition of the appropriate amount of peptide. Incubate at room temperature for 1 hr. Centrifuge both tubes at 10,000Xg for 5 min. Use the supernatants from tubes 1 and 2 for parallel experiments.

**STORAGE:** Lyophilized antibody and lyophilized peptide can be stored intact at room temperature for several weeks. For longer periods they should be stored at -20°C. After reconstitution, the antibody solution can be stored at 4°C for up to 2 weeks. For longer periods, small aliquots should be stored at -20°C or below. Avoid freeze/thaw cycles. Further dilutions should be made using a carrier protein such as 1% BSA. Centrifuge all antibody preparations before use (10,000 X g for 5 minutes). After reconstitution, control peptide should be stored at -20°C.

## REFERENCES:

1. A. Moroni *et al.* *J Biol Chem.* 2001 **276** 29233
2. W. Han *et al.* *Circ Res.* 2002 **91** 790
3. T.N. Doan *et al.* *J Neurosci.* 2004 **24** 3335
4. V. Nacri and A.A. Accili. *J Biol Chem.* 2004 **279** 16832
5. M. Jiang *et al.* *Circulation.* 2004 **109** 1783
6. B. Much *et al.* *J Biol Chem.* 2003 **278** 43781

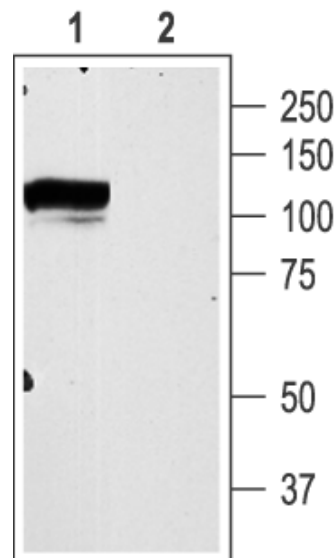


Figure: Lanes 1 & 2: Rat brain membranes. Lane 1 was probed with Anti-HCN2. Lane 2 was probed with Anti-HCN2 which had been pre-incubated with the control peptide.

FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.