



6405 Mira Mesa Blvd, Suite 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

## **Data Sheet**

### **CNGA2 Expression Vector** **Catalog #: 60642**

**Description:** The CNGA2 expression vector is designed to express human cyclic nucleotide-gated cation channel 2 (CNGA2) in mammalian cells.

**Application:** Suitable for use in transient assays with the PDE1B-HEK293 Cell Line (BPS Bioscience, #60402).

**Size:** 20 µg

**Concentration:** 80 ng/µl

**Storage buffer:** 10 mM Tris-HCl (pH 8), 1 mM EDTA

**Storage condition:** Store plasmid at -20°C in aliquots. Avoid repeated freeze/thaw cycles.

**Stability:** Stable for at least one year when stored as recommended.

**Note:** The vector is suitable for transient transfection. It is not suitable for amplification in bacteria.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**  
Or you can Email us at: [\*\*support@bpsbioscience.com\*\*](mailto:support@bpsbioscience.com)  
Please visit our website at: [\*\*www.bpsbioscience.com\*\*](http://www.bpsbioscience.com)