



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

ANPRA/Aequorin Expression Vector Catalog #: 60641

Description: The ANPRA/Aequorin expression vector is designed to co-express human atrial natriuretic peptide receptor A (ANPRA, also called natriuretic peptide receptor A/guanylate cyclase A) and jellyfish (*Aequorea victoria*) Aequorin in mammalian cells.

Application: Suitable for use with the PDE1B Cell-based Reporter Assay Kit, Cat. # 60601.

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 **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com