

# Di-2-ANEPEQ [JPW 1114]

Catalog number: 21496

Unit size: 5 mg

#### **Product Details**

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

#### **Chemical Properties**

Appearance Solid

Molecular Weight 549.38

Soluble In Water

Chemical Structure

### **Spectral Properties**

Excitation Wavelength 488 nm

Emission Wavelength 701 nm

## **Applications**

Di-2-ANEPEQ is used for monitoring fast membrane potential changes. ANEP dyes belong to the class of the fast-response membrane potential dyes. Their optical response is fast enough to detect transient membrane potential changes in excitable cells where they demonstrate a membrane potential-dependent shift in excitation spectra. This feature allows the measurement of membrane potential changes by excitation ratio. These dyes are weakly fluorescent in aqueous media, and become strongly fluorescent upon binding to lipophilic environments (such as membranes). In general, fast-response probes operate by means of a change in their electronic structure, and consequently their fluorescence properties, in response to a change in the surrounding electric field. Their optical response is sufficiently fast to detect transient (millisecond) potential changes in excitable cells, including single neurons, cardiac cells and intact brains. However, the magnitude of their potential-dependent fluorescence change is often small; fast-response probes typically show a 2-10% fluorescence change per 100 mV.