



Product Data Sheet

Product Name:	Oxonol VI [<i>Bis - (3 - propyl - 5 - oxoisoxazol - 4 - yl)pentamethine oxonol</i>]
Size:	100 mg
Catalog Number:	AS-84704
Molecular Weight:	316.4
Solvent:	DMSO
Product Description:	Oxonol VI is used as an optical indicator for membrane potentials in lipid vesicles. Oxonol VI is a sensitive slow-response membrane potential probe whose magnitude of optical response is much larger than that of faster response probes (Oxonol V). The fluorescence of Oxonol VI decreases upon membrane hyperpolarization. In general, slow response probes are suitable for detecting changes in average membrane potentials of nonexcitable cells caused by respiratory activity, ion-channel permeability, drug binding and other factors.
Spectra:	Maximum Ex/Em wavelength = 599/634 nm
Shelf Life and Storage:	Shelf life is two years if stored at -20C, desiccated and protected from light.
Product citation	Mennerick, S. et al. (2010). Diverse Voltage-Sensitive Dyes Modulate GABAAR Receptor Function. <i>J Neurosci</i> 30, 2871.
Related Products:	

This product is for *in vitro* research use only