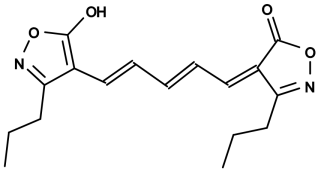
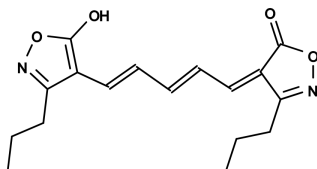


**Oxonol VI [Bis-(3-propyl-5-oxoisoxazol-4-yl)
pentamethine oxonol]**Catalog number: 21417
Unit size: 25 mg**Product Details**

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	316.35
Soluble In	DMSO
Chemical Structure	

**Applications**

Oxonol VI is a sensitive slow-response membrane potential probe that has faster response than Oxonol V. The fluorescence of Oxonol VI decreases upon membrane hyperpolarization. In general, slow-response probes exhibit potential-dependent changes in their transmembrane distribution that are accompanied by a fluorescence change. The magnitude of their optical responses is much larger than that of fast-response probes (typically a 1% fluorescence change per mV). Slow-response probes, which include cationic carbocyanines, rhodamines and anionic oxonols, are suitable for detecting changes in average membrane potentials of nonexcitable cells caused by respiratory activity, ion-channel permeability, drug binding and other factors.