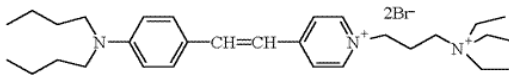
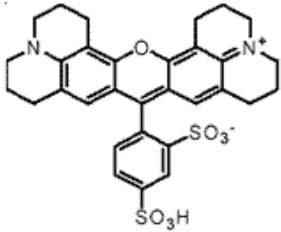


Instruction Manual

| | |
|--------------------------|--|
| Catalog Number | PK-CA707-70032 |
| Description | The Nerve Terminal Green and Red Staining Kits have been designed to detect and study recycling of the synaptic vesicles in neuronal synapses and neuromuscular junctions. Although each kit uses a different amphiphilic styryl pyridinium dye to visualize the vesicles, preparation of the tissue, the required reagents and the staining procedures are identical for each. The dyes included in the kits fluoresce intensely inside plasma membranes but only minimally in aqueous environments. When neurons are actively releasing neurotransmitters, dye is incorporated into recycled vesicles at the presynaptic terminal. The dye-containing vesicles are not susceptible to subsequent wash steps that remove the excess dye from the outer leaflet of the plasma membrane. Thus, only active presynaptic terminals are labeled. |
| Quantity | 1 set (5x1 mg SynaptoGreen C4 and 100 mg Sulforhodamine 101) |
| SynapseGreen C4 | |
| Description | A widely used green fluorescent dye for following synaptic activities by staining synaptic vesicles at the synapse or neuromuscular junctions. ¹ When used in combination with the red fluorescent dye SynapseRed PK-CA707-70021), synapses or neuromuscular junctions can be imaged independently in two colors. Recent research results from Tsien's ² and Kay's ³ labs showed that brain slices can be stained with the styryl dye SynapseGreen when SR101 (Sulforhodamine 101, PK-CA707-80101) or ADVASEP-7 (PK-CA707-70029) is used to reduce the background staining. PromoKine offers several Nerve Terminal Staining Kits for staining brain slices. |
| Molecular Structure |  |
| Molecular Formula | C ₃₀ H ₄₉ Br ₂ N ₃ |
| Molecular Weight | 612 g/mol |
| Appearance | Red solid |
| Purity | ≥95% (determined by HPLC) |
| Absorption/Emission max. | λ _{ex} λ _{em} : 510/625 nm (in MeOH); 480/598 nm (in membranes). |
| Solubility | Soluble in water |
| References | <ol style="list-style-type: none"> 1. J. Neurosci. 12, 363(1992) 2. Science. 255, 200(1992) 3. Neuron 24, 803(1999) 4. Neuron 24, 809(1999) |
| Intended Use | For in vitro research use only. Not for diagnostic or therapeutic procedures. |
| Storage | Store desiccated at 4°C. |

| | |
|---------------------------------|---|
| Sulforhodamine 101 | |
| Description | Sulforhodamine 101 has a longer absorption wavelength than sulforhodamine B. Like sulforhodamine B and G, these fluorescent dyes have been primarily used as polar tracers for the studies of neuronal cell morphology and cell-cell communications - in addition to their potential use in cancer drug screening. |
| Molecular Structure |  <p>The chemical structure of Sulforhodamine 101 consists of a central xanthenoquinone core. The core is substituted with two piperazine rings, one on each side of the oxygen atom in the xanthene moiety. The nitrogen atom in the left piperazine ring is neutral, while the nitrogen atom in the right piperazine ring is positively charged. The central carbon of the xanthene ring is substituted with a 4,4'-disulfonatophenylene group, which consists of a benzene ring with two sulfonate groups (-SO₃⁻ and -SO₃H) at the para positions.</p> |
| Molecular Formula | C ₃₁ H ₃₀ N ₂ O ₇ S ₂ |
| Molecular Weight | 607 g/mol |
| Appearance | Red solid |
| Absorption/Emission max. | λ _{ex} \ λ _{em} : 586/605 nm |
| Solubility | Soluble in water |
| Intended Use | For in vitro research use only. Not for diagnostic or therapeutic procedures. |
| Storage | Store desiccated at RT. |

PromoCell GmbH

Sickingenstr. 63/65
69126 Heidelberg
Germany

Email: info@promokine.info
www.promokine.info

North America

Phone: 1 - 866 - 251 - 2860 (toll free)
Fax: 1 - 866 - 827 - 9219 (toll free)

Deutschland

Telefon: 0800 - 776 66 23 (gebührenfrei)
Fax: 0800 - 100 83 06 (gebührenfrei)

France

Téléphone: 0800 90 93 32 (ligne verte)
Téléfax: 0800 90 27 36 (ligne verte)

United Kingdom

Phone: 0800 - 96 03 33 (toll free)
Fax: 0800 - 169 85 54 (toll free)

Other Countries

Phone: +49 6221 - 649 34 0
Fax: +49 6221 - 649 34 40