Revised: June 05, 2018

Product Information

SynaptoGreen™ & SynaptoRed™ Nerve Terminal Dyes

| Terriman byes | | | | | | |
|-------------------|--|----------|----|----|--------|--|
| Catalog Number | Product | Size | m* | n* | MW | |
| 70042 | - SynaptoGreen™ C1 | 5 mg | 0 | 1 | 527.4 | |
| 70043 | | 5 x 1 mg | | | | |
| 70044 | SynaptoGreen™ C2 (equivalent to FM®2-10) | 5 mg | 1 | 1 | 556 | |
| 70045 | | 5 x 1 mg | | | | |
| 70023 | - SynaptoGreen™ C3 | 5 mg | 2 | 1 | 542 | |
| 70026 | | 5 x 1 mg | | | | |
| 70020 | SynaptoGreen™ C4 (equivalent to FM®1-43) | 5 mg | 3 | 1 | 612 | |
| 70022 | | 5 x 1 mg | | | | |
| 70046 | SynaptoGreen™ C5 (equivalent to FM®1-84) | 5 mg | 4 | 1 | 639.59 | |
| 70047 | | 5 x 1 mg | | | | |
| 70048 | SynaptoGreen™ C18 (equivalent to FM®3-25) | 5 mg | 17 | 1 | 1004 | |
| 70049 | | 5 x 1 mg | | | | |
| 70040 | - SynaptoRed™ C1 | 5 mg | 0 | 3 | 580 | |
| 70041 | | 5 x 1 mg | | | | |
| 70021 | SynaptoRed™ C2 (equivalent to FM®4-64) | 5 mg | 1 | 3 | 608 | |
| 70027 | | 5 x 1 mg | | | | |
| 70019 | SynaptoRed™C2M (equivalent to FM®5-95) | 5 x 1 mg | 1 | 3 | 565.43 | |
| 70028 | | 5 mg | | | | |
| | | | | | | |

^{*} See Figure 1.

Storage and Handling

Store desiccated at 4°C or below. Protect from light, especially in solution. Product is stable for at least 12 months from date of receipt when stored as recommended. SynaptoGreen and SynaptoRed dyes are soluble in water. Stock solutions can be prepared at 10 mM and stored at 4°C or -20°C for six months or longer.

Spectral Properties

SynaptoGreen dyes:

Abs/Em (MeOH): 510/625 nm

Abs/Em (membranes): ~480/600 nm (Fig. 2)

SynaptoRed dyes:

Abs/Em (MeOH): 543/- nm (emission in MeOH is too weak to measure) Abs/Em (membranes): ~510/750 nm (Fig. 3)

Note: Excitation/emission settings of 515/640 nm have been used for detection of yeast vacuole staining with SynaptoRed C2 (FM4-64) (Ref. 1); these settings should be applicable to SynaptoRed C1.

$$[{\rm CH_3}({\rm CH_2})_{\rm m}]_2{\rm N} - ({\rm CH=CH})_{\rm n} - ({\rm CH_2})_3 - {\rm N}({\rm CH_2CH_3})_3$$

Figure 1. General structure of SynaptoGreen and SynaptoRed dyes, where m = 0-17 and n = 1 or 3. The positively charged head group of SynaptoRed C2M is a trimethylammonium group instead of a triethylammonium group.

Product Description

Nerve terminal probes are a series of fluorescent cationic styryl dyes developed to follow synaptic activities at neuromuscular junctions or synapses. These dyes typically have a lipophilic tail (two carbon chains) at one end and a highly hydrophilic, cationically charged head group at the other end. These nerve terminal probes were originally called FM® dyes, and are available from Biotium under the trademark names of SynaptoGreen $^{\text{TM}}$ and SynaptoRed $^{\text{TM}}$. SynaptoGreen probes are dyes with a single double bond (n = 1) while SynaptoRed probes are dyes with three double bonds (n = 3). A nerve terminal dye is named as either SynaptoGreen or SynaptoRed followed by a carbon number designating the length of the lipophilic tail

Cationic styryl dyes are believed to function by staining synaptic vesicles in an activity-dependent fashion. In the presence of cells or tissue preparations, the dyes partition between the aqueous phase, where the dyes are virtually non-fluorescent, and the outer leaflet of the cell surface membranes, where the dyes insert the lipophilic end into the membranes and become intensely fluorescent. During endocytosis following nerve stimulation, the dyes become trapped inside the vesicles. Thus, after washing off the dyes on the cell surface, the fluorescent signal is proportional to the number of newly formed vesicles. On the other hand, during exocytosis, the dyes are released from the vesicles along with neurotransmitters, causing a decrease in fluorescent signal. As a result, the change in fluorescent intensity reflects the amount of endocytosis/exocytosis or synaptic activity. The rate of fluorescence increase during exocytosis, the "on-rate", and the rate of fluorescence decrease during exocytosis, the "off-rate", vary from dye to dye. In general, dyes with longer lipophilic tails and more double bonds have a higher affinity toward membrane and thus a higher on-rate and lower off-rate.

Some styryl dyes can enter cells through ion channels. SynaptoGreen C18, which has a long carbon chain and cannot pass through ion channels, has been used as a control to distinguish mechanisms of dye uptake (2).

SynaptoGreen and SynaptoRed are not fixable. The AM and HM dyes (see Related Products) are aldehyde-fixable nerve terminal dyes.

When using nerve terminal dyes, one frequent problem researchers encounter is background fluorescence due to nonspecific membrane staining. Although most of the background fluorescence can be removed by repeated washing, the problem is still significant with dyes that have a longer tail or more double bonds, particularly when the dyes are used in tissue preparations. Biotium offers various background reducing agents for use with nerve terminal dyes. ADVASEP-7 is a sulfonated beta-cyclodextrin that helps to remove dye during washing (3). SCAS quenches fluorescence of membrane-bound dye without the need for repeated wsh steps (4). Sulforhodamine 101 is a red fluorescent dye that quenches background fluorescence from green nerve terminal dyes by FRET (5). We also offer Nerve Terminal Staining Kits that include dyes and background reducing agents (see Related Products).

References

- 1. Vida, TA and Emr, SD. J. Cell Biol 128, 779(1995).
- 2. Meyers, JR, et al. J Neurosci., 23, 4054(2003).
- 3. Kay, AR, et al. Neuron 24, 809 (1999).
- 4. Höltje, M, et al. J. Biol. Chem. 283(14):9289(2008)
- 5. Pyle, JL, et al. Neuron 24, 803 (1999).

Continued next page

Assay Protocol

The following is an example of a protocol for nerve terminal staining of cultured neurons on coverslips. Nerve terminal dyes also can be used to label endocytic vesicles in non-neuronal cell types. Staining can be performed at 4°C for selective labeling of the plasma membrane; at room temperature or 37°C, endocytosis of the dye generally occurs within 10 minutes. Buffers other than Tyrode solution may be used. The addition of the sodium channel blocker tetrodotoxin (TTX) is optional, its purpose is to block action potentials and prevent synaptic vesicle release after staining. Optimal protocols for specific applications may need to be determined by the user; see references for examples of protocols for staining brain slices and other cell types.

- Dilute nerve terminal dye to a final concentration of 4 uM in 50 mM Tyrode solution. Place the coverslip with your cells in this solution for 1 minute at room temperature. Use enough solution to completely submerge the cells.
- Transfer the coverslip to Tyrode + 0.5 uM tetrodotoxin (TTX, catalog no. 00061) solution for 1 minute at room temperature.
- 3. Wash the coverslip several times in Tyrode + TTX at room temperature.
 Note: to reduce background, 1 mM ADVASEP-7 (catalog no. 70029) can be added to the wash solution. Alternatively, SCAS (catalog no. 70037) can be used to quench background without repeated washes. Incubate the coverslip for 4 minutes at room temperature in Tyrode + TTX + 0.5 mM SCAS.
- 4. Mount the coverslip in Tyrode + TTX and image.

Note: for SynaptoGreen dyes, 50 uM sulforhodamine 101 (catalog no. 80101) can be included during mounting to quench extracellular fluorescence.

Note: SynaptoGreen and SynaptoRed dyes are not fixable. The AM and HM dyes (see Related Products) are aldehyde-fixable nerve terminal dyes.

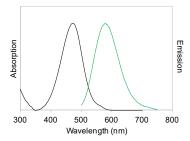


Figure 2. Absorption and emission spectra of SynaptoGreen C4 in liposomes. Spectra for other SynaptoGreen dyes are similar.

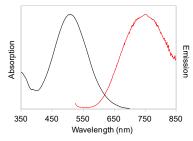


Figure 3. Absorption and emission spectra of SynaptoRed C2 in liposomes. Spectra for other SynaptoRed dyes are similar.

Related Products

| Catalog number | Product | | |
|----------------|--|--|--|
| 70024 | AM1-43 | | |
| 70038 | AM1-44 | | |
| 70036 | AM2-10 | | |
| 70051 | AM3-25 | | |
| 70025 | AM4-64 | | |
| 70039 | AM4-65 | | |
| 70050 | AM4-66 | | |
| 70029 | ADVASEP-7 | | |
| 70037 | SCAS | | |
| 80101 | Sulforhodamine 101 | | |
| 70030 | Nerve Terminal Staining Kit I 5 x 1 mg SynaptoGreen™ C4 and 250 mg ADVASEP-7 | | |
| 70031 | Nerve Terminal Staining Kit II (A) 1 mg AM1-43 and 100 mg ADVASEP-7 | | |
| 70031-1 | Nerve Terminal Staining Kit II (B) 1 mg AM1-43 and 100 mg SCAS | | |
| 70032 | Nerve Terminal Staining Kit III 5 x 1 mg SynaptoGreen™ C4 and 100 mg Sulforhodamine 101 | | |
| 70034 | Nerve Terminal Staining Kit V 5 x 1 mg SynaptoRed™ C2 and 250 mg ADVASEP-7 | | |
| 00060 | Tetrodotoxin, citrate-free | | |
| 00061 | Tetrodotoxin, with citrate | | |
| 00010 | α -Bungarotoxin | | |
| 00019 | β-Bungarotoxin | | |

Please visit our website at www.biotium.com for information on our life science research products, including fluorescent CF™ dye bungarotoxins, antibodies, and other conjugates, calcium and other ion indicator dyes, apoptosis detection reagents, and other fluorescent probes and kits for cell biology research.

FM is a registered trademark of Thermo Fisher Scientific.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.