

ADVASEP-7

ADVASEP-7 is a sulfonated β -cyclodextrin derivative that has been reported to reduce background fluorescence when using SynptoGreen C4 (also called FM1-43, catalog no. [70020](#)) to stain brain slices.



Product attributes

Product Description

ADVASEP-7 is a sulfonated beta-cyclodextrin derivative that has been reported to reduce background fluorescence when using SynptoGreen™ C4 (also called FM@1-43, catalog no. [70020](#)) to stain brain slices.

- Average MW: 2163
- $C_{42}H_{70-n}O_{35}(C_4H_8SO_3Na)_n \cdot n$ hydrate where n = Average Degree of Substitution 6.5
- White solid soluble in water
- Store at 4 °C

Nerve Terminal Dyes and Kits

Product	Catalog No.	Features
SynptoGreen™ Dyes (Ex/Em ~480/660 nm in membranes)	70020... 70053	View full list of SynptoGreen™ Dyes
SynptoRed Dyes™ (Ex/Em ~510/750 nm in membranes)	70021... 70050	View full list of SynptoRed™ Dyes
Background Reducers		
ADVASEP-7	70029	Sulfonated cyclodextrin that aids in removal of free dye during washes
SCAS	70037	Quenches extracellular fluorescences with fewer wash steps than ADVASEP-7
Sulforhodamine 101	80101	Red fluorescent dye that quenches extracellular fluorescence of SynptoGreen™ dyes
Nerve Terminal Staining Kits		
Nerve Terminal Staining Kit I	70030	Includes SynptoGreen™ C4 and ADVASEP-7
Nerve Terminal Staining Kit II (A)	70031	Includes AM1-43 and ADVASEP-7
Nerve Terminal Staining Kit II (B)	70031-1	Includes AM1-43 and SCAS
Nerve Terminal Staining Kit III	70032	Includes SynptoGreen™ C4 and Sulforhodamine 101
Nerve Terminal Staining Kit V	70034	Includes SynptoRed™ C2 and ADVASEP-7

m is the number of carbons in the lipophilic tail and n is the number of double bonds linking the two aromatic rings in the dye.

**The positively-charged end of SynptoRed C2M is a trimethylammonium group.FM is a registered trademark of Thermo Fisher Scientific.

References

1. Neuron 24, 809 (1999).

This datasheet was generated on January 7, 2026 at 04:37:38 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/advasep-7/>