

# Product Information

## Hydroxystilbamidine

Product	Catalog No.	Unit Size
Hydroxystilbamidine, methanesulfonate	80014	10 mg
Hydroxystilbamidine, 4% in H <sub>2</sub> O	80023	200 uL

### Storage and Handling

Store at 4°C, protected from light. Product is stable for at least 12 months from date of receipt when stored as recommended.

### Molecular Information

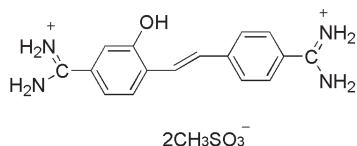
C<sub>18</sub>H<sub>24</sub>N<sub>4</sub>O<sub>7</sub>S<sub>2</sub>

MW: 473

CAS: 223769-64-0

Color & Form: Yellow solid (80014); Yellow liquid (80023)

Solubility: Soluble in water at 4% (solubility in PBS buffer is significantly lower)



### Spectral Properties

Absorption: ~290-370 nm (broad absorption in water)

Ex/Em: When bound to membranes at pH 7.4, excitation: 350-395 nm (broad); emission: 530-600 nm (broad).

### Product Description

Hydroxystilbamidine (or Fluoro-Gold™) has been used extensively as a retrograde tracer for neurons and also a histochemical stain. The dye has very broad spectra with absorption from 350-400 nm and emission from 530-620 nm. A wide band UV excitation filter is recommended for optimal excitation.

Hydroxystilbamidine, 4% solution in deionized H<sub>2</sub>O is a convenient, ready-made solution. If desired, the solution can be diluted to a lower concentration. For most of the applications, the typical dye concentration is 2% to 4%. Note: The dye solution is produced using type I ultrapure water, but it is not prepared or packaged aseptically, and should be sterile-filtered for aseptic applications. We offer Mini-Syringe Filters and Mini-Syringe Filtration Kit with syringes and sterile vials (see Related Products), which are convenient for 0.2 um filtration of small volumes of aqueous solutions.

### References

1) Brain Res. 377, 147(1986); 2) Brain Res. 475, 259(1988); 3) J. Neurosci. Methods, 23, 1(1988); 4) Brain Res. 553, 135(1991); 5) J. Neurosci. Methods, 45, 87(1992).

### Related Products

Catalog number	Product
22025	Mini Syringe Filters (Set of 5)
22026	Mini Syringe Filtration Kit (for 5 Samples)
70020	SynaptoGreen™ C4 (equivalent to FM@1-43)
70021	SynaptoRed™ C2 (equivalent to FM@4-64)
80027	PathoGreen™ Histofluorescent Stain
80028	Congo Red, High Purity Grade
80030	DCDAPH
80033	Thioflavin T, High Purity Grade
00060	Tetrodotoxin, Citrate-Free
00061	Tetrodotoxin, With Citrate
00010	α-Bungarotoxin
23001	EverBrite™ Mounting Medium
23002	EverBrite™ Mounting Medium with DAPI
23003	EverBrite™ Hardset Mounting Medium
23004	EverBrite™ Hardset Mounting Medium with DAPI
23005	CoverGrip™ Coverslip Sealant
23007	TrueBlack™ Lipofuscin Autofluorescence Quencher
23012	TrueBlack® IF Background Suppressor System (Permeabilizing)
40061	RedDot™2 Far Red Nuclear Counterstain
22005	Mini Super <sup>HT</sup> Pap Pen 2.5 mm tip, ~400 uses
22006	Super <sup>HT</sup> Pap Pen 4 mm tip, ~800 uses
22020	10X Phosphate-Buffered Saline (PBS)
22023	Paraformaldehyde, 4% in PBS, Ready-To-Use Fixative

Please visit our website at [www.biotium.com](http://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, neuronal tracers, apoptosis detection reagents, and other fluorescent probes and kits for cell biology research.

Fluoro-Gold is a trademark of Fluorochrome, LLC.

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