

Hydroxystilbamidine (Fluoro-Gold™), 4% in H₂O

Hydroxystilbamidine (also called Fluoro-Gold™) has been used extensively as a retrograde tracer for neurons and also a histochemical stain. The 4% solution in deionized H₂O is a convenient, ready-made solution.



Product attributes

CAS number	223769-64-0
Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Green
Excitation/Emission	361/536 nm

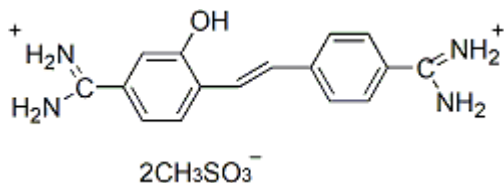
Product Description

Hydroxystilbamidine (also called Fluoro-Gold™) has been used extensively as a retrograde tracer for neurons and also a histochemical stain. Fluoro-Gold™ can undergo retrograde axonal transport, it shows extensive filling of dendrites, and has a high resistance to fading.

This product is provided as a ready-made 4% solution in deionized H₂O. If desired, the solution can be diluted to a lower concentration. For most applications, the typically 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](#), which are convenient for 0.2 um filtration of small volumes of aqueous solutions.

We also sell hydroxystilbamidine in solid form (10 mg, cat#[80014](#)).

- $\lambda_{Ex}/\lambda_{Em} = 361/536$ nm
- Store at 4 °C and protect from light
- C₁₈H₂₄N₄O₇S₂
- MWt: 473
- [223769-64-0]



Fluoro-Gold is a trademark of Fluorochrome, LLC.

References

1. Brain Res 553, 135 (1991).
2. Brain Res 377, 147 (1986)