Glowing products for science

## 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 H<sub>2</sub>O 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

Call us: 800-304-5357 Email: btinfo@biotium.com

## **Product Description**

Hydroxystilbamidine (also called Fluoro-Gold<sup>TM</sup>) has been used extensively as a retrograde tracer for neurons and also a histochemical stain. Fluoro-Gold<sup>TM</sup> 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  $H_2O$ . 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).

- $\lambda Ex/\lambda Em = 361/536 \text{ nm}$
- Store at 4°C and protect from light
- C<sub>18</sub>H<sub>24</sub>N<sub>4</sub>O<sub>7</sub>S<sub>2</sub>
- MWt: 473
- [223769-64-0]

2CH<sub>3</sub>SO<sub>3</sub>

Fluoro-Gold is a trademark of Fluorochrome, LLC.

## References

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

•

This datasheet was generated on November 6, 2025 at 12:33:18 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/hydroxystilbamidine-4-in-h2o-equivalent-to-fluorogoldtm/">https://biotium.com/product/hydroxystilbamidine-4-in-h2o-equivalent-to-fluorogoldtm/</a>