

Hydroxystilbamidine (Fluoro-Gold™)

Hydroxystilbamidine (also called Fluoro-Gold $^{\text{TM}}$) has been used extensively as a retrograde tracer for neurons and also a histochemical stain.



Product attributes

Call us: 800-304-5357

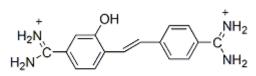
CAS number	223769-64-0
Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Green
Excitation/Emission	361/536 nm

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.

Also see hydroxystilbamidine 4% in H_2O (80023), which is a more convenient, ready-to-use form.

- $\lambda Ex/\lambda Em = 361/536 \text{ nm}$
- Yellow solid soluble in water
- Store at 4°C and protect from light
- C₁₈H₂₄N₄O₇S₂
- MW: 473
- [223769-64-0]



2CH3SO3

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 23, 2025 at 10:44:55 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/hydroxystilbamidine-equivalent-to-fluorogoldtm/