

Hydroxystilbamidine (Fluoro-Gold™)

Hydroxystilbamidine (also called Fluoro-Gold™) has been used extensively as a retrograde tracer for neurons and also a histochemical stain.



Product attributes

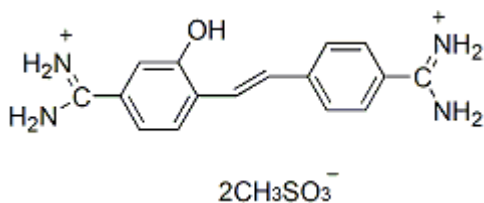
CAS number	223769-64-0
Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Green
Excitation/Emission	361/536 nm

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.

Also see hydroxystilbamidine 4% in H₂O ([80023](#)), which is a more convenient, ready-to-use form.

- λEx/λEm = 361/536 nm
- Yellow solid soluble in water
- Store at 4 °C and protect from light
- C₁₈H₂₄N₄O₇S₂
- MW: 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).

This datasheet was generated on June 21, 2026 at 12:15:27 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/hydroxystilbamidine-equivalent-to-fluorogoldtm/>