

# Hydroxystilbamidine (ultra pure)

## Neuronal detection

Hydroxystilbamidine is a cationic dye, equivalent to Fluoro- Gold™, which is frequently used as a retrograde neuronal tracer. The fluorescent tracer provides intense retrograde labeling that is extremely sensitive and reliable, does not diffuse out of retrograde-labeled neurons, and can be pressure-injected or introduced by iontophoresis into cells. Hydroxystilbamidine is compatible with most frequently used neuroanatomical techniques such as immunofluorescence, immunocytochemistry, autoradiography, and horseradish peroxidase-based histochemistry, paraffin embedding. Wavelength Maxima: Excitation 385nm, Emission 536nm

Citations: 4

[View Online »](#)

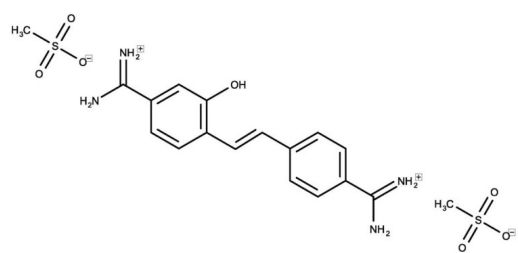
## Ordering Information

[Order Online »](#)

ENZ-52253	10mg
-----------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored as recommended.
Handling	Protect from light. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	4-[(E)-2-(4-Carbamimidoylphenyl)ethenyl]-3-hydroxybenzenecarboximidamide
CAS	223769-64-0 (methanesulfonate salt)
Formula	$C_{18}H_{24}N_4O_7S_2$
MW	472.5
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.

Technical Info / Product Notes	This product is a member of the CELLESTIAL <sup>®</sup> product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. CELLESTIAL <sup>®</sup> reagents and kits are optimal for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required.
--------------------------------	---



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)