



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

PathoGreen™ Histofluorescent Stain, 1000X in water

PathoGreen™ Histofluorescent Stain is an anionic green fluorescent dye functionally similar to Fluoro-Jade dyes. These dyes stain degenerating neurons and their processes after exposure to a variety of neurotoxic insults in brain sections and cultured neurons.



Product attributes

Apoptosis/viability marker	Dead cell stain
For live or fixed cells	For fixed cells/tissue sections
Detection method/readout	Fluorescence microscopy
Assay type/options	Endpoint assay
Colors	Green
Excitation/Emission	497/520 nm

Product Description

PathoGreen™ Histofluorescent Stain is an anionic green fluorescent dye (Ex/Em 497/520 nm) functionally similar to Fluoro-Jade® dyes. These dyes stain degenerating neurons and their processes after exposure to a variety of neurotoxic insults in brain sections and cultured neurons. The mechanism of neuronal staining by anionic fluorescent dyes has not been determined. It has been proposed that the negatively charged dyes bind to positively charged polyamines or other molecules specifically generated in dying neurons.

Fluoro-Jade is a registered trademark of Histo-Chem, Inc.

References

Shimada, R. et al. (2014). Neurological Research 36(4), 239-246.
[doi:10.1179/1743132813Y.0000000289](https://doi.org/10.1179/1743132813Y.0000000289)

This datasheet was generated on June 10, 2026 at 09:10:18 PM. Visit [product page](https://biotium.com/product/pathogreen-histofluorescent-stain-1000x-in-water/) to check for updated information before use.
Product link: <https://biotium.com/product/pathogreen-histofluorescent-stain-1000x-in-water/>