Lucifer Yellow Cadaverine Biotin-X, dipotassium salt

This fluorescent biotin derivative is a useful tracer for studying neuronal morphology. The biotin portion of the molecule allows researchers to amplify the fluorescent signal using standard biotin-avidin or biotin-streptavidin biochemistry.



Product attributes

Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Green
Excitation/Emission	428/532 nm
Conjugation	Biotin

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

Product Description

This fluorescent biotin derivative is a useful tracer for studying neuronal morphology. The biotin portion of the molecule allows researchers to amplify the fluorescent signal using standard biotin-avidin or biotin-streptavidin biochemistry. The staining can tolerate aldehyde fixation, sectioning, and washing with detergent.

- λ_{Abs}/λ_{Ems} 428/532 nm
- Yellow solid soluble in water
- Store desiccated at ≤4 °C and protect from light
- C33H42K2N6O11S3
- MW: 873

References

- 1. J Neurosci Meth 53, 23 (1994).
- 2. J Neurosci Meth 46, 59 (1993).

This datasheet was generated on November 12, 2025 at 01:40:15 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/lucifer-yellow-cadaverine-biotin-x-dipotassium-salt/