

## 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.



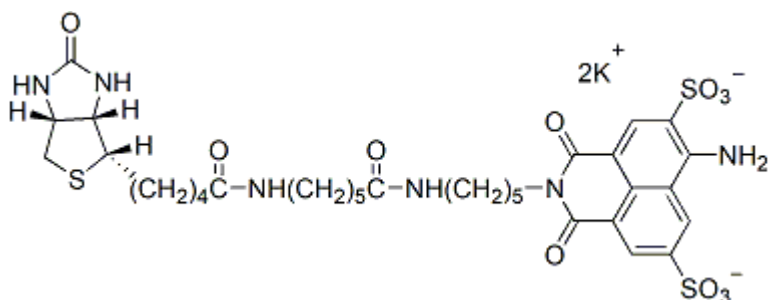
### Product attributes

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

### 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.

- $\lambda_{Abs}/\lambda_{Ems}$  428/532 nm
- Yellow solid soluble in water
- Store desiccated at  $\leq 4^{\circ}\text{C}$  and protect from light
- $\text{C}_{33}\text{H}_{42}\text{K}_2\text{N}_6\text{O}_{11}\text{S}_3$
- MW: 873



### References

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

This datasheet was generated on June 22, 2026 at 04:59:00 AM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/lucifer-yellow-cadaverine-biotin-x-dipotassium-salt/>