## Lucifer Yellow CH, lithium salt, 100 mM in water

Biotum

Lucifer Yellow CH lithium salt (LY CH lithium salt) is a widely used polar molecular tracer for studying neuronal morphology. The fluorescent molecule contains a carbohydrazide that allow the molecule to be aldehyde-fixable.

## CAS number

Product attributes	
CAS number	67769-47-5
Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Green
Excitation/Emission	428/536 nm

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

## **Product Description**

Lucifer Yellow CH lithium salt (LY CH lithium salt) is a widely used polar molecular tracer for studying neuronal morphology. The fluorescent molecule contains a carbohydrazide that allow the molecule to be aldehyde-fixable. The lithium salt form of Lucifer Yellow is in an aqueous solution format that is ready for microinjection.

- $\lambda_{Ex}/\lambda_{Em}=428/536$  nm
- Yellow solution. Store at room temperature and protect from light.
- C<sub>13</sub>H<sub>9</sub>Li<sub>2</sub>N<sub>5</sub>O<sub>9</sub>S<sub>2</sub>
- MW: 457
- [67769-47-5]

This datasheet was generated on November 11, 2025 at 05:09:25 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/lucifer-yellow-ch-lithium-salt-100mm-in-water/">https://biotium.com/product/lucifer-yellow-ch-lithium-salt-100mm-in-water/</a>