## Lucifer Yellow CH, lithium salt

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.



## **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.

- $\lambda_{Ex}/\lambda_{Em} = 428/536 \text{ nm}$
- Yellow solid soluble in water
- Store at 4°C 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]

## References

- 1. Nature 292, 17 (1981).
- 2. Cell 14, 741 (1978).

This datasheet was generated on November 17, 2025 at 11:08:25 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/lucifer-yellow-ch-lithium-salt/">https://biotium.com/product/lucifer-yellow-ch-lithium-salt/</a>