

Lucifer Yellow CH, lithium salt, 100 mM in water

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.



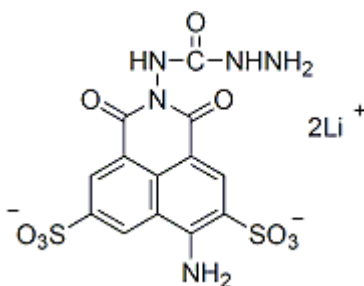
Product attributes

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

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.
- $C_{13}H_9Li_2N_5O_9S_2$
- MW: 457
- [67769-47-5]



This datasheet was generated on June 9, 2026 at 08:02:45 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/lucifer-yellow-ch-lithium-salt-100mm-in-water/>