Lucifer Yellow CH, potassium salt

Lucifer Yellow CH potassium salt (LY CH potassium 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	71206-95-6
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 potassium salt (LY CH potassium 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.

- $\bullet \ \lambda_{Ex}/\lambda_{Em} = 428/536 \ nm$
- Yellow solid soluble in water
- Store at 4°C and protect from light
- \bullet C₁₃H₉K₂N₅O₉S₂
- MW: 522
- [71206-95-6]

References

1. Nature 292, 17 (1981).

2. Cell 14, 741 (1978).

This datasheet was generated on November 11, 2025 at 05:09:24 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/lucifer-yellow-ch-potassium-salt/