Lucifer Yellow Cadaverine

Like Lucifer Yellow CH, Lucifer Yellow Cadaverine can be used as a fixable fluorescent tracer. In addition, the dye can be used to conjugate to carboxylic acid groups.



Product attributes

CAS number	149733-79-9
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

Like Lucifer Yellow CH, Lucifer Yellow Cadaverine can be used as a fixable fluorescent tracer. In addition, the dye can be used to conjugate to carboxylic acid groups.

- λ_{Ex}/λ_{Em}= 428/536 nm
- Yellow solid soluble in water
- Store at 4°C and protect from light
- C₁₇H₁₇K₂N₃O₈S₂
- MW: 534

References

1. Anal Biochem 211, 210 (1993).

This datasheet was generated on November 8, 2025 at 01:38:51 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/lucifer-yellow-cadaverine/