Biocytin

Biocytin (ϵ -Biotinoyl-L-lysine) is a very useful cellular tracer that can be introduced into cells by microinjection.



Product attributes

CAS number	576-19-2
Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Conjugation	Biotin

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

Product Description

Biocytin (ϵ -Biotinoyl-L-lysine) is a very useful cellular tracer that can be introduced into cells by microinjection.

- White solid soluble in water or DMSO
- Store at 4°C
- C₁₆H₂₈N₄O₄S
- MW: 372.48
- [576-19-2]

References

- 1. Anat Embryol 199, 149 (1998).
- 2. Brain Res 782, 113 (1998).
- 3. Brain Res Devel Brain Res 106, 79 (1998).
- 4. Exp Brain Res 119, 375 (1998).

This datasheet was generated on November 11, 2025 at 06:55:37 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/biocytin-e-biotinoyl-l-lysine/