

## Biocytin

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



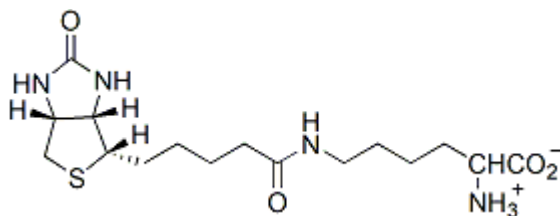
### Product attributes

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

## 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
- C<sub>16</sub>H<sub>28</sub>N<sub>4</sub>O<sub>4</sub>S
- 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 January 12, 2026 at 06:03:04 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/biocytin-e-biotinoyl-l-lysine/>