

## Biocytin hydrazide

Biocytin hydrazide is an aldehyde-fixable cellular tracer that can be introduced into cells by microinjection. Biocytin hydrazide can also be used to label glycoproteins or other carbohydrate molecules (See [Reactive Dyes for Labeling Glycoproteins and Other Carbohydrate Molecules](#)).



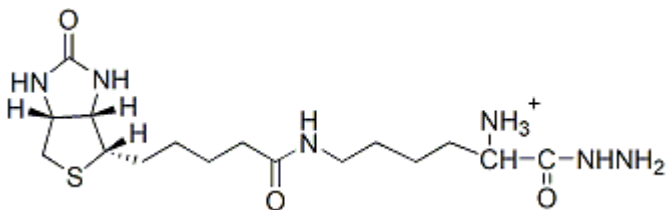
### Product attributes

CAS number	102743-85-1
------------	-------------

## Product Description

Biocytin hydrazide is an aldehyde-fixable cellular tracer that can be introduced into cells by microinjection. Biocytin hydrazide can also be used to label glycoproteins or other carbohydrate molecules (See [Reactive Dyes for Labeling Glycoproteins and Other Carbohydrate Molecules](#)).

- White solid soluble in DMSO
- Store desiccated at 4 °C
- C<sub>16</sub>H<sub>30</sub>N<sub>6</sub>O<sub>3</sub>S
- MW: 386.51
- [102743-85-1]



## References

1. Glycoconjugate J 9, 21 (1992).
2. Anal Biochem 171, 1 (1988).
3. Anal Biochem 170, 271 (1988).
4. Biochem Biophys Res Comm 136, 80 (1986).

This datasheet was generated on June 3, 2026 at 07:02:32 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/biocytin-hydrazide/>