## **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

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

## **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
- $\bullet$  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 November 5, 2025 at 02:59:13 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/biocytin-hydrazide/">https://biotium.com/product/biocytin-hydrazide/</a>