

## Fluorescein biotin

Fluorescein biotin (5-((N-(5-(N-(6-(Biotinoyl)amino)hexanoyl)amino)pentyl)thioureidyl) fluorescein) can be used to detect and quantify biotin binding sites and the degree of biotinylation of proteins. Also see biotin-rhodamine 110 ([80022](#)) and biotin-4-fluorescein ([90062](#)).



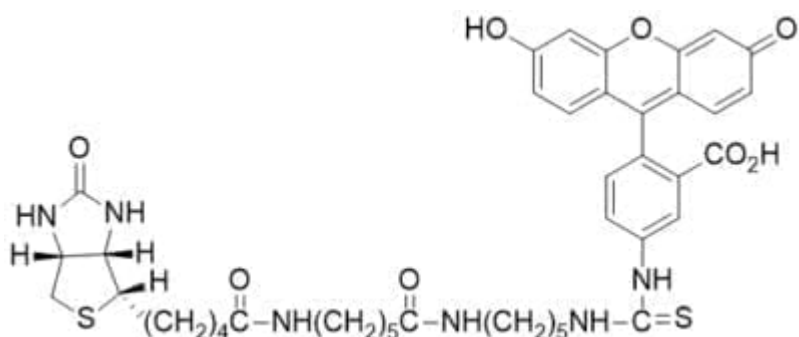
### Product attributes

|                             |                     |
|-----------------------------|---------------------|
| CAS number                  | 134759-22-1         |
| Probe cellular localization | Fluid phase tracer  |
| Cell permeability           | Membrane impermeant |
| Colors                      | Green               |
| Excitation/Emission         | 494/518 nm          |
| Conjugation                 | Biotin              |

## Product Description

Fluorescein biotin (5-((N-(5-(N-(6-(Biotinoyl)amino)hexanoyl)amino)pentyl)thioureidyl) fluorescein) can be used to detect and quantify biotin binding sites and the degree of biotinylation of proteins. Also see biotin-rhodamine 110 ([80022](#)) and biotin-4-fluorescein ([90062](#)).

- $\lambda_{Ex}/\lambda_{Em}$  = 494/518 nm
- Yellow solid soluble in DMSO, or DMF
- Store at -20 °C and protect from light
- $C_{42}H_{50}N_6O_8S_2$
- MW: 831
- [134759-22-1]



## References

1. Biochim Biophys Acta 1381, 203 (1998).
2. Clin Chem 40, 2112 (1994).

This datasheet was generated on December 23, 2025 at 03:15:01 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/fluorescein-biotin-5-n-5-n-6-biotinoylaminohexanoylaminopentylthioureidyl-fluorescein/>