

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](#)).



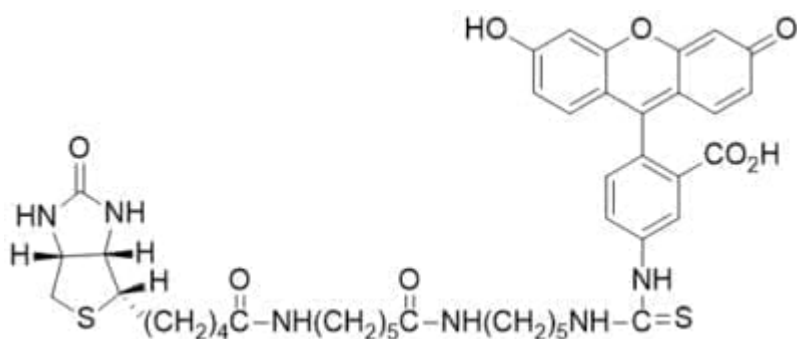
Product attributes

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
- C₄₂H₅₀N₆O₈S₂
- MW: 831
- [134759-22-1]



References

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

This datasheet was generated on June 18, 2026 at 09:02:24 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/>