

# Product Information

## Biotin-4-Fluorescein

**Catalog Number:** 90062

**Unit Size:** 10 mg

### Storage and Handling

Store desiccated at -20°C. Protect from light. Product is stable for at least 12 months from date of receipt when stored as recommended.

**Molecular Information:** C<sub>33</sub>H<sub>32</sub>N<sub>4</sub>O<sub>8</sub>S

**CAS number:** 1032732-74-3

**Molecular Weight:** 645

**Color and Form:** Yellow solid

**Purity:** >95%

**Solubility:** Soluble in DMSO, or pH>7 buffers

**Absorption/Emission:** 494/523 nm

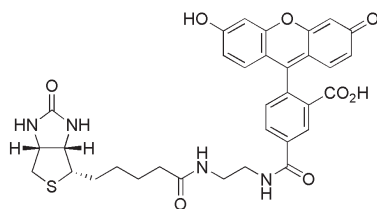


Figure 1. Biotin-4-fluorescein

### Product Description

Biotin-4-fluorescein has been reported to be a superior fluorescent biotin derivative for accurate measurement of avidin and streptavidin in crude biological samples. Binding of biotin to avidin and streptavidin is high affinity, fast and noncooperative.

[Browse our website](#) to view our full selection of biotin and CF® Dye conjugates.

### References

- PNAS 102, 24, pp 8507 (2005);
- Biochim Biophys Acta 1427, 33 (1999);
- Biochim Biophys Acta 1427, 44 (1999).

### Related Products

Cat. No.	Product
92208...92228	CF® Dye & Biotin SE Protein Labeling Kits
00017	Biotin-XX-α-Bungarotoxin
00020	Biotin-cAMP
60022	Biotin-DHPE
80019	Fluorescein Biotin
80022	Biotin-Rhodamine 110
92167	Biotin Azide
92176	Biotin-XX Tyramide
92168	Biotin Alkyne
29030... 29086	CF® Dye Streptavidin Conjugates
29049	HRP Streptavidin
29071	Alkaline Phosphatase Streptavidin
29048	APC Streptavidin
29044	R-PE Streptavidin
40061	RedDot™2 Far Red Nuclear Counterstain, 200X in DMSO
40083...41040	NucSpot® Nuclear Stains
23001	EverBrite™ Mounting Medium
23002	EverBrite™ Mounting Medium with DAPI
23003	EverBrite™ Hardset Mounting Medium
23004	EverBrite™ Hardset Mounting Medium with DAPI
23005	CoverGrip™ Coverslip Sealant
22015	Fixation Buffer, 100 mL
22016	Permeabilization Buffer, 100 mL
22017	Permeabilization and Blocking Buffer, 100 mL
22010	10% Fish Gelatin Blocking Buffer, 100 mL
22011	Fish Gelatin Powder, 2 X 50 g
22014	30% Bovine Serum Albumin Solution, 100 mL
22002	Tween®-20, 50 mL

Please visit our website at [www.biotium.com](http://www.biotium.com) for information on our life science research products, including fluorescent CF® Dye antibody conjugates and reactive dyes, biotinylation reagents with a full selection of reactive functional groups, as well as easy-to-use protein and antibody biotinylation kits, and kits for cell biology research.

CF dye technology is covered by pending US and international patents. TWEEN is a registered trademark of CRODA AMERICAS LLC.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.