

## Flubi-2

Flubi-2 is the hydrolyzed product of the site-directed fluorogenic probe Flubida-2 ([#51022](#)) and can be used as a standard for calibration.



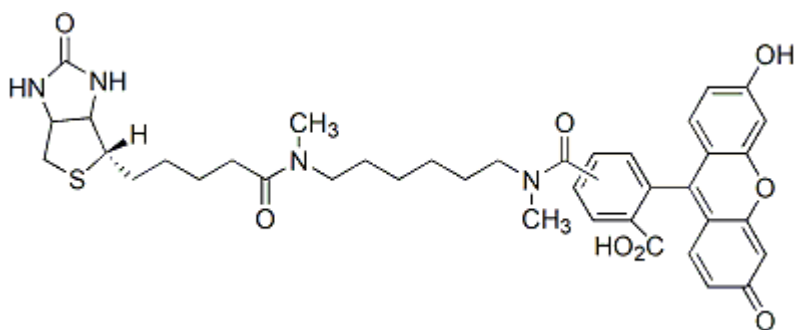
### Product attributes

Cell permeability	Membrane impermeant
pKa	pKa 6.5
Indicator type	Ratiometric
Colors	Green
Excitation/Emission	492/517 nm

## Product Description

Flubi-2 is the hydrolyzed product of the site-directed fluorogenic probe Flubida-2 ([#51022](#)) and can be used as a standard for calibration. The dye can also be microinjected into cells to probe the local pH where avidin fusion proteins are located.

- $\lambda_{Ex}/\lambda_{Em}$ : 492/517 nm (pH 9)
- Yellow solid soluble in DMSO
- Store at -20 °C
- $C_{39}H_{44}N_4O_8S$
- MW: 728.85



## References

1. Chemistry and Biology 7, 197 (2000).

This datasheet was generated on January 7, 2026 at 03:23:31 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/flubi-2/>