## 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
рКа	pKa 6.5
Indicator type	Ratiometric
Colors	Green
Excitation/Emission	492/517 nm

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

## **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.

- λ<sub>Ex</sub>/λ<sub>Em:</sub> 492/517 nm (pH 9)
- Yellow solid soluble in DMSO
- Store at -20°C
- C<sub>39</sub>H<sub>44</sub>N<sub>4</sub>O<sub>8</sub>S
- MW: 728.85

## References

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

This datasheet was generated on November 2, 2025 at 04:56:01 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/flubi-2/">https://biotium.com/product/flubi-2/</a>