

Glowing products for science

H2DCFDA

H₂DCFDA (2',7'-Dichlorodihydrofluorescein diacetate) is a useful fluorogenic reagent to detect reactive oxygen intermediates in cells. On oxidation, H₂DCFDA becomes the highly green fluorescent 2',7'-dichlorofluorescein.



Call us: 800-304-5357 Product attributes

CAS number 4091-99-0

Excitation/Emission 504/529 nm (end product)

Product Description

 H_2DCFDA (2',7'-Dichlorodihydrofluorescein diacetate) is a useful fluorogenic reagent to detect reactive oxygen intermediates in cells. On oxidation, H_2DCFDA becomes the highly green fluorescent 2',7'-dichlorofluorescein.

- $\lambda_{Ex}/\lambda_{Em}$ (pH >6) = 504/529 nm (end product)
- White solid soluble in DMSO or DMF
- Store desiccated at -20°C and protect from light
- C₂₄H₁₆Cl₂O₇
- MW: 487.29
- [4091-99-0]

References

- 1. Plant Growth Regul (2010) doi: 10.1007/s10725-010-9545-y
- 2. Physiologia Plantarum (2010) doi: 10.1111/j.1399-3054.2010.01400.x
- 3. Biochemistry 39, 1040 (2000).
- 4. Cancer Res. 60, 219 (2000).
- 5. Free Rad Res 32, 57 (2000).
- 6. J Immunol Meth 159, 173 (1993).
- 7. J Immunol Meth159, 131 (1993).
- 8. Cytometry 13, 615 (1992).

This datasheet was generated on November 26, 2025 at 04:48:33 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/27-dichlorodihydrofluorescein-diacetate-h2dcfda/