

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.



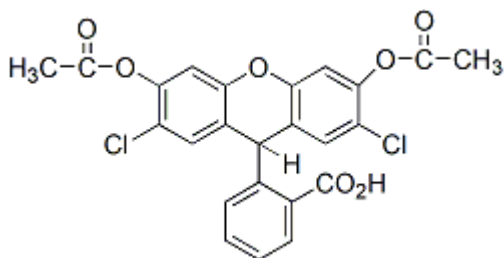
Product attributes

CAS number	4091-99-0
Excitation/Emission	504/529 nm (end product)

Product Description

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.

- $\lambda_{Ex}/\lambda_{Em}$ (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 Meth 159, 131 (1993).
8. Cytometry 13, 615 (1992).

This datasheet was generated on June 14, 2026 at 10:31:36 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/27-dichlorodihydrofluorescein-diacetate-h2dcfda/>