

DCDAPH (DANIR-2c)

Far-red fluorescent probe with high affinity for A β plaques/aggregates, that can be used for fluorescent staining of brain sections and *in vivo* detection of A β ₁₋₄₂ aggregates in small animals.



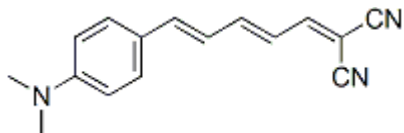
Product attributes

Apoptosis/viability marker	Amyloid stain
Colors	Far-red, Near-infrared
Excitation/Emission	597/665 nm

Product Description

DCDAPH (1,1-dicyano-6-(4-N,N-dimethylaminophenyl)-1,3,5-hexatriene) aka DANIR-2c, can be used for fluorescent staining of brain sections, as well as *in vivo* detection of A β ₁₋₄₂ aggregates by small animal near-IR imaging¹.

- Far-red fluorescent probe with high affinity for A β plaques/aggregates
- Affinity for A β ₁₋₄₂ aggregates: K_i= 37 nM, K_d=27 nM
- Dark purple solid
- $\lambda_{Ex}/\lambda_{Em}$ = 597/665 nm (PBS)
- Store at -20 °C and protect from light, especially in solution
- C₁₆H₁₅N₃
- MW: 249.31



References

1. J. Phys. Chem. B DOI: 10.1021/acs.jpcc.2c07783
2. J. Am. Chem. Soc. dx.doi.org/10.1021/ja4052922

This datasheet was generated on January 25, 2026 at 01:40:56 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/dcdaph/>