

RH421

RH421 is a fast responding membrane potential dye, exhibiting >20% fluorescence change per 100 mV on neuroblastoma cells.



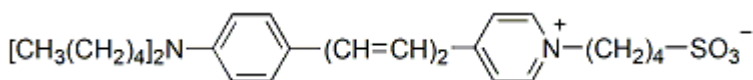
Product attributes

CAS number	107610-19-5
For live or fixed cells	For live/intact cells
Potential dependence	Fast-response membrane potential dye
Colors	Far-red
Excitation/Emission	515/704 nm

Product Description

RH421 is a fast responding membrane potential dye, exhibiting >20% fluorescence change per 100 mV on neuroblastoma cells.

- $\lambda_{Ex}\lambda_{Em}$ = 515/704 nm (MeOH)
- Red solid soluble in DMSO, EtOH or DMF
- Store at -20 °C and protect from light, especially in solution
- $C_{29}H_{42}N_2O_3S$
- MW: 498.72
- [107610-19-5]



References

1. Biophys J 42, 195 (1983).

This datasheet was generated on December 21, 2025 at 04:04:06 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/rh421/>