RH421

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



Product attributes

| . Todaot attributor | |
|-------------------------|--------------------------------------|
| 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 |

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

Product Description

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

- λεχλεm= 515/704 nm (MeOH)
- Red solid soluble in DMSO, EtOH or DMF
- Store at -20°C and protect from light, especially in solution
- C₂₉H₄₂N₂O₃S
- MW: 498.72
- [107610-19-5]

$$[CH_3(CH_2)_4]_2N$$
 $-(CH=CH)_2$ N $-(CH_2)_4$ $-SO_3$

References

1. Biophys J 42, 195 (1983).

This datasheet was generated on November 2, 2025 at 05:08:12 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/rh421/