RuBP-2S

RuBP-2S

(Ruthenium(bis(bathophenanthroline))(bathophenanthrolinedisulfonate) is a luminescent ruthenium complex similar to RuBPS (80021).



Product attributes

Probe cellular localization	Fluid phase tracer
Cell permeability	Membrane impermeant
Colors	Red
Excitation/Emission	462 (broad)/603 nm

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

Product Description

RuBP-2S (Ruthenium(bis(bathophenanthroline))(bathophenanthrolinedisulfonate) is a luminescent ruthenium complex similar to RuBPS (80021). Two of the bathophenanthroline ligands are neutral while the third bathophenanthroline ligand is bis-sulfonated, resulting in an overall neutral charge of the ruthenium complex. RuBP-2S shares the spectral properties of RuBPS and may also have similar applications.

- $\lambda_{Ex}/\lambda_{Em}(H2O) = 462 \text{ (broad)/603 nm}$
- Yellow solid soluble in water
- Store at 4°C and protect from light
- C72H46N6O6RuS2
- MW: 1256.3

This datasheet was generated on November 1, 2025 at 11:20:28 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/rubp-2s-rutheniumbisbathophenanthrolinedisulfonate/