

RuBP-2S

RuBP-2S (Ruthenium(bis(bathophenanthroline))(bathophenanthroline disulfonate)) 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

Product Description

RuBP-2S (Ruthenium(bis(bathophenanthroline))(bathophenanthroline disulfonate)) 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.

- $\lambda_{Ex}/\lambda_{Em}(H_2O) = 462 \text{ (broad)}/603 \text{ nm}$
- Yellow solid soluble in water
- Store at 4 °C and protect from light
- $C_{72}H_{46}N_6O_6RuS_2$
- MW: 1256.3

