

## RuBP-4S

RuBP-4S ([Ruthenium(bathophenanthroline)(bis(bathophenanthrolinedisulfonate))],disodium salt) 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-4S ([Ruthenium(bathophenanthroline)(bis(bathophenanthrolinedisulfonate))],disodium salt) is a luminescent ruthenium complex similar to RuBPS ([80021](#)). Two of the bathophenanthroline ligands are bis-sulfonated while the third bathophenanthroline ligand is neutral, resulting in an overall double negative charges of the ruthenium complex. RuBP-4S shares the spectral properties of RuBPS and may also have similar applications.

- $\lambda_{Ex}/\lambda_{Em}(H_2O) = 462$  (broad)/603 nm
- Yellow solid soluble in water
- Store at 4 °C and protect from light
- $C_{72}H_{44}N_6Na_2O_{12}RuS_4$
- MW: 1460.46

