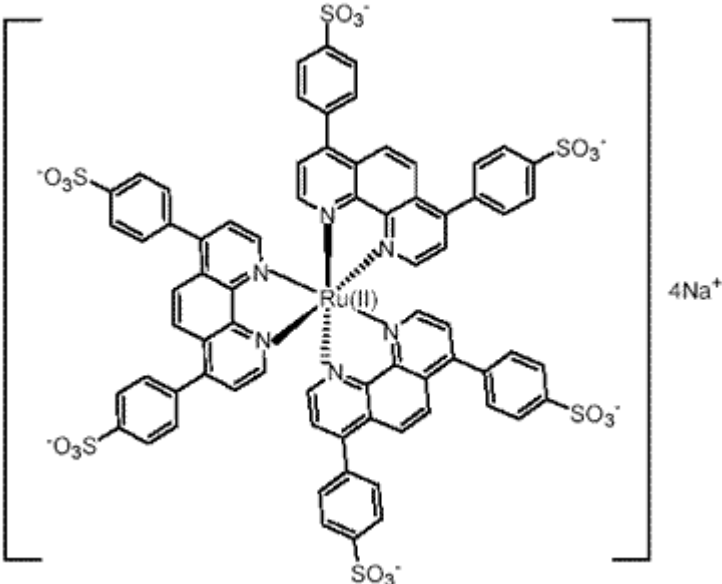


## Ruthenium(tris(bathophenanthrolinedisulfonate), tetrasodium salt

## Instruction Manual

Catalog Number	PK-CA707-80021
Description	RuBPS is a negatively-charged water-soluble luminescent ruthenium complex. The complex can be efficiently excited by either UV light or visible light in the wavelength range from 400 nm to 500 nm with bright red emission centered around 605 nm. Ruthenium complexes such as RuBPS typically have luminescence lifetimes in the micro second range and a Stokes shift over 100 nm. By using time-resolved luminescence measurement, the combination of long luminescent lifetime and large Stokes shift offer the possibility for extremely sensitive detection that could approach the sensitivity of radioactive detection. Among many potential applications, RuBPS could be used as a polar tracer similar to the uses of Lucifer Yellow and FluoroGold.
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex}$ \ $\lambda_{em}$ : 462 (broad)/603 (in water) nm; Extinction coefficient = 32,000
Molecular Structure	
Molecular Weight / Molecular Formula	1664.55 Da; C <sub>72</sub> H <sub>42</sub> N <sub>6</sub> Na <sub>4</sub> O <sub>18</sub> RuS <sub>6</sub>
Purity	NA
Appearance / Formulation / Solubility	Yellow solid; soluble in water.
Storage & Stability	Store at 4°C. Protect from light.
Applications	See Description
References	NA
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.