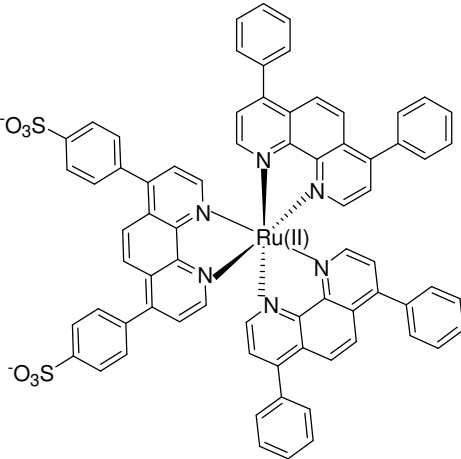


(Ruthenium(bis(bathophenanthroline))(bathophenanthroline disulfonate))

Instruction Manual

Catalog Number	PK-CA707-80024
Description	RuBP-2S is a luminescent ruthenium complex similar to RuBPS (Cat.No. PK-CA707-80021). 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 (see Cat.No. PK-CA707-80021).
Quantity	5 mg
Excitation / Emission Maxima	$\lambda_{ex} \setminus \lambda_{em}$: 462 (broad)/603 (in MeOH) nm; Extinction coefficient = 16,000
Molecular Structure	
Molecular Weight / Molecular Formula	1256.3 Da; C ₇₂ H ₄₆ N ₆ O ₆ RuS ₂
Purity	NA
Appearance / Formulation / Solubility	Yellow solid; soluble in water and DMF.
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.