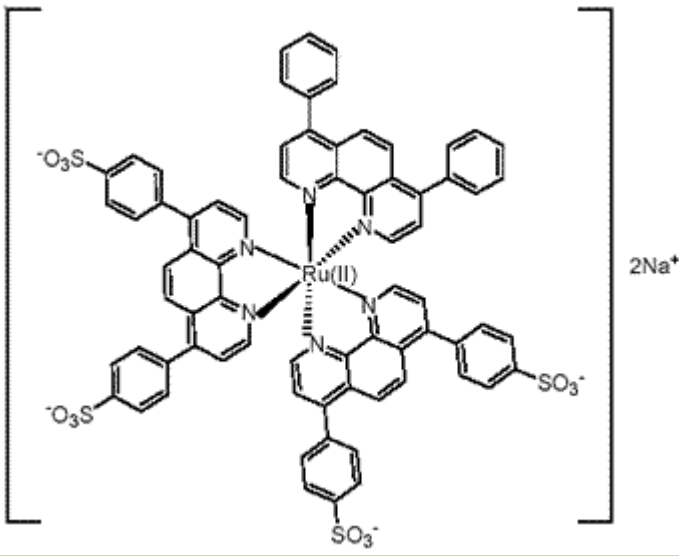


[Ruthenium(bathophenanthroline)(bis(bathophenanthrolinedisulfonate))], disodium salt**Instruction Manual**

| | |
|---------------------------------------|--|
| Catalog Number | PK-CA707-80025 |
| Description | RuBP-4S is a luminescent ruthenium complex similar to RuBPS (Cat.No. PK-CA707-80021). 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 (see also Cat.No. PK-CA707-80021). |
| Quantity | 5 mg |
| Excitation / Emission Maxima | $\lambda_{ex} \setminus \lambda_{em}$: 462 (broad)/603 (in water) nm; Extinction coefficient = 32,000 |
| Molecular Structure |  |
| Molecular Weight / Molecular Formula | 1460.46 Da; $C_{72}H_{44}N_6Na_2O_{12}RuS_4$ |
| 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.