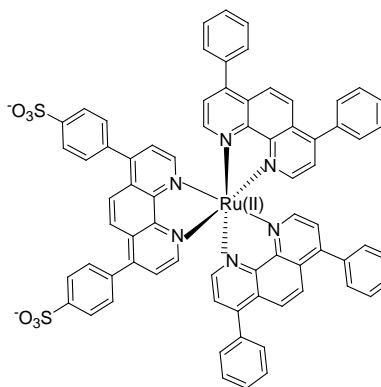


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** **RuBP-2S** (Ruthenium(bis(bathophenanthroline))(bathophenanthroline disulfonate))**CATALOG #** **80024****MOLECULAR INFORMATION:**  $C_{72}H_{46}N_6O_6RuS_2$   
MWt: 1256.3**PROPERTIES:**

**Color & Form** Yellow solid  
**Solubility** Soluble in DMF  
**Absorption/Emission** 462 (broad)/603 nm (in MeOH)  
**Extinction Coefficient** 16,000

**STORAGE AND HANDLING:** Store at 4 °C and protect from light**APPLICATION:** RuBP-2S 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 (See product description for #80021).**RELATED PRODUCTS** *RuPBS (#80021); RuBP-4S(#80025)***TOXICITY:** Not established. Not listed by NTP, IARC or OSHA.

<b>FIRST AID:</b>	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.
-------------------	---

<b>Disclaimer:</b> <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--