

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

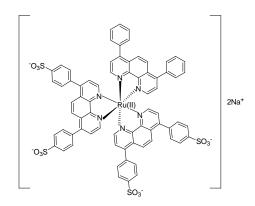
## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** RuBP-4S ([Ruthenium(bathophenanthroline)(bis(bathophenanthrolinedisulfonate))],

disodium salt)

**CATALOG #** 80025

**INFORMATION:** MWt: 1460.46



PROPERTIES:

**Color & Form** Yellow solid **Solubility** Soluble in water

**Absorption/Emission** 462 (broad)/603 nm (in H<sub>2</sub>O)

Extinction 32,000

Coefficient

STORAGE AND Store at 4 °C and protect from light

HANDLING:

**APPLICATION:** RuBP-4S 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 (See product description for #80021).

**TOXICITY:** Not established. Not listed by NTP, IARC or OSHA.

FIRST AID: 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.

**Disclaimer:** 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.