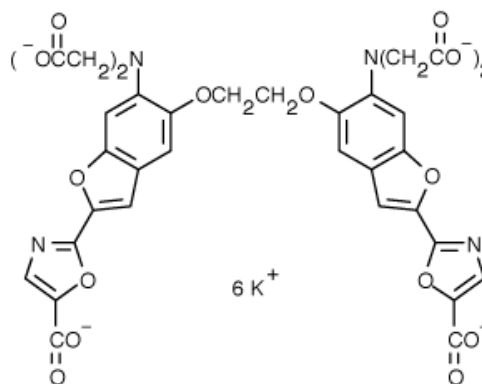


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Bis-Fura-2, hexapotassium salt**CATALOG #:** 50045**MOLECULAR INFORMATION:** C<sub>34</sub>H<sub>20</sub>K<sub>6</sub>N<sub>4</sub>O<sub>18</sub>  
MWt: 1007.14**PROPERTIES:**

<b>Color &amp; Form</b>	Light yellow solid
<b>Purity</b>	≥ 98% by TLC
<b>Solubility</b>	Soluble in water (pH >6)
<b>Absorption/Emission</b>	363 nm/512 nm (no Ca <sup>2+</sup> ); 335 nm/505 nm (high Ca <sup>2+</sup> )
<b>Extinction Coefficient</b>	56,000 M <sup>-1</sup> cm <sup>-1</sup> (363nm, no Ca <sup>2+</sup> ); 68,000 M <sup>-1</sup> cm <sup>-1</sup> (high Ca <sup>2+</sup> )

**STORAGE AND HANDLING:**

Store desiccated at 4°C or -20 °C. Protect from light. It can be reconstituted in aqueous buffers having pH >6 and stored at 4°C. Aqueous solution has a shelf life of at least 3 months if stored at 4°C or below.

**APPLICATION:**

Bis-Fura-2, hexapotassium salt is membrane-impermeant derivative of membrane-permeant calcium indicator Fura-2, AM ester. It contains two Fura-2-pype fluorophores. This indicator has a slightly lower infinity for Ca<sup>2+</sup> and can be introduced into cell by microinjection.

Ref: 1) *J Neurophysiol* **84**, 2777(2000); 2) *Brain Res* **831**, 113(1999); 3) *J Neurophysiol* **81**, 2508(1999); 4) *J Neurosci Res* **57**, 906(1999); 5) *J. Neurochem Int* **34**, 391(1999)

**TOXICITY:**

Unknown

**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.