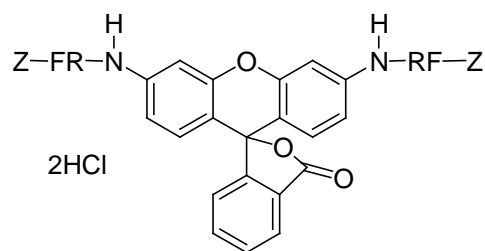


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Rhodamine 110, bis-(*N*-CBZ-*L*-phenylalanyl-*L*-arginine amide), dihydrochloride (Z-FR-R110)

**CATALOG #:** 10209

**MOLECULAR INFORMATION:** C<sub>66</sub>H<sub>72</sub>Cl<sub>2</sub>N<sub>12</sub>O<sub>11</sub>  
MW: 1278.26

**PROPERTIES:**

**Color & Form** Pinkish red solid  
**Purity** ≥ 75% by HPLC  
**Solubility** Soluble in DMSO  
**Absorption/Emission** 496 nm/520 nm (strong fluorescence)  
(end product R110)

**STORAGE AND HANDLING:** Store desiccated at ≤ 4 °C. Protect from light.

**APPLICATION:** Z-FR-R110 is a substrate for the cysteine proteases cathepsins B and L. The substrate has been used in flow cytometry studies of human monocytes and rat macrophages.

Ref.: 1) *Glia* **7**, 183(1993); 2) *Biol Chem Hoppe Seyler* **373**, 433(1992); 3) *Biol Chem Hoppe Seyler* **373**, 547(1992)

**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.