



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Z-FR-R110

This item is discontinued.



Product attributes

Excitation/Emission	496/520 nm (end product)
---------------------	--------------------------

Product Description

This item is discontinued.

References

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

This datasheet was generated on January 17, 2026 at 11:15:22 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/z-fr-r110-rhodamine-110-bis-n-cbz-l-phenylalanyl-l-arginine-amide-dihydrochloride/>