BZiPAR

BZiPAR (Rhodamine 110,

bis-(N-CBZ-L-isoleucyl-L-prolyl-L-arginine amide), dihydrochloride) is a substrate for trypsin (1,2). The substrate has been reported to enter live cells and be hydrolyzed by lysosomal proteases (3).

sfortum #80

Product attributes

Excitation/Emission

496/520 nm (end product)

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

BZiPAR (Rhodamine 110, bis-(N-CBZ-L-isoleucyl-L-prolyl-L-arginine amide), dihydrochloride) is a substrate for trypsin (1,2). The substrate has been reported to enter live cells and be hydrolyzed by lysosomal proteases (3).

- $\lambda_{Ex}\lambda_{Em}$ of end product (R110) = 496/520 nm
- Off-white to pink solid soluble in DMSO
- Store at 4°C and protect from light
- C₇₀H₈₈Cl₂N₁₄O₁₃
- MW: 1404.5

References

- 1. PNAS 97, 13126 (2000).
- 2. Biochim Biophys Acta 788, 74 (1984).
- 3. Photochem Photobiol 44, 461 (1986).

This datasheet was generated on November 2, 2025 at 05:53:12 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/bzipar-rhodamine-110-bis-n-cbz-l-isoleucyl-l-prolyl-l-arginine-amide-dihydrochloride/