

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).



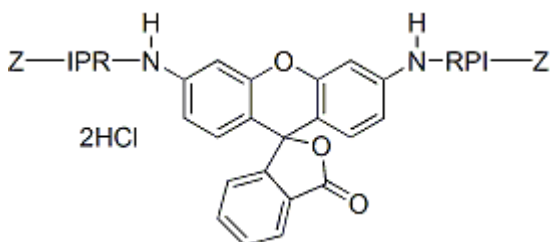
Product attributes

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

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
- $C_{70}H_{88}Cl_2N_{14}O_{13}$
- 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 June 13, 2026 at 06:09:47 PM. 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/>