

BZAR

BZAR (Rhodamine 110, bis-(N-CBZ-L-arginine amide), dihydrochloride) is a substrate for serine proteases (1-5) and is also an inhibitor for quatinidinobenzoatase. It is 50 to 100 times more sensitive than the AMC-based substrate.



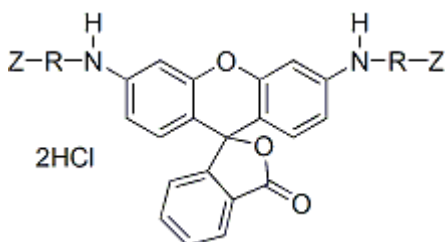
Product attributes

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

Product Description

BZAR (Rhodamine 110, bis-(N-CBZ-L-arginine amide), dihydrochloride) is a substrate for serine proteases (1-5) and is also an inhibitor for quatinidinobenzoatase. It is 50 to 100 times more sensitive than the AMC-based substrate.

- $\lambda_{Ex}\lambda_{Em}$ of end product (R110) = 496/520 nm
- Off-white to pink solid soluble in DMSO
- Store at 4 °C and protect from light
- $C_{48}H_{52}Cl_2N_{10}O_9$
- MW: 983.91



References

1. J Histochem Cytochem 49, 1473(2001).
2. Biotechniques 29, 1108-1113 (2000).
3. J Enzyme Inhib 6, 303(1992).
4. Biochem Soc Trans 18, 632 (1990).
5. J Cancer Res Clin Oncol 116, 57 (1990).
6. J Enzyme Inhib 2, 209 (1988).

This datasheet was generated on December 6, 2025 at 11:22:59 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/bzar-rhodamine-110-bis-n-cbz-l-arginine-amide-dihydrochloride/>