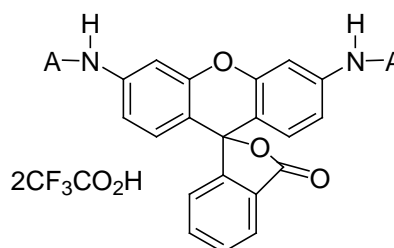


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Rhodamine 110, bis-(L-aspartic acid amide), dinitrfluoroacetic acid salt (D₂R)

CATALOG #: 10210

MOLECULAR INFORMATION: C₃₂H₂₆F₆N₄O₁₃
MW: 788.57

**PROPERTIES:**

Color & Form Off-white to pinkish solid
Purity ≥ 95% by TLC
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: D₂R is a general substrate for caspases including caspases 3 and 7. Unlike other fluorogenic substrates for caspases, D₂R appears to be membrane-permeant, and thus can be used to measure caspase activity in live cells without osmotic shock for flow cytometry studies, as well as for cell extracts.

Ref.: 1) *Biochemistry* **38**, 13906(1999).

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: <i>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.</i>
--