

D2R

D₂R (Rhodamine 110, bis-(L-aspartic acid amide), ditrifluoroacetic acid salt) 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.



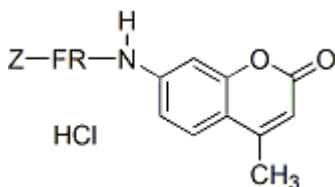
Product attributes

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

Product Description

D₂R (Rhodamine 110, bis-(L-aspartic acid amide), ditrifluoroacetic acid salt) 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. For real-time detection of caspase-3 activity in intact cells, see our novel [NucView@488 Caspase-3 Substrate](#).

- Ex/Em of end product (R110) = 496/520 nm
- Off-white to pink solid soluble in DMSO
- Store at 4 °C and protect from light
- C₃₂H₂₆F₆N₄O₁₃
- MW: 788.57



References

1. Biochemistry 38, 13906 (1999).

This datasheet was generated on December 6, 2025 at 12:24:50 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/d2r-rhodamine-110-bis-l-aspartic-acid-amide-ditrifluoroacetic-acid-salt/>