

PRODUCT INFORMATION



(D)₂-Rh 110 (trifluoroacetate salt) Item No. 28213

Formal Name: (3S,3'S)-4,4'-[(3-oxospiro[isobenzofuran-1(3H),9'-[9H]xanthene]-3',6'-diyl)diimino]bis[3-amino-4-oxo-butanoic acid, trifluoroacetate salt

Synonyms: D₂R, (Asp)₂-Rhodamine 110, Rhodamine 110 bis-(L-aspartic acid amide)

MF: C₂₈H₂₄N₄O₉ • XCF₃COOH

FW: 560.5

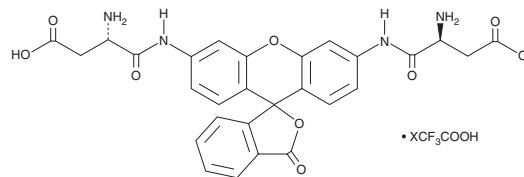
Purity: ≥95%

Ex./Em. Max: 496/520 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

(D)₂-Rh 110 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the (D)₂-Rh 110 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. (D)₂-Rh 110 (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

(D)₂-Rh 110 is a fluorogenic caspase substrate.¹ Upon enzymatic cleavage by caspases, rhodamine 110 (Item No. 19061) is released in an apoptosis-dependent manner and its fluorescence can be used to quantify caspase activity. Rhodamine 110 displays excitation/emission maxima of 496/520 nm, respectively.

Reference

1. Hug, H., Los, M., Hirt, W., et al. Rhodamine 110-linked amino acids and peptides as substrates to measure caspase activity upon apoptosis induction in intact cells. *Biochemistry* **38(42)**, 13906-13911 (1999).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/05/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM