PRODUCT INFORMATION



Rhodamine 110 (chloride)

Item No. 19061

CAS Registry No.:	13558-31-1	\sim
Formal Name:	3,6-diamino-9-(2-carboxyphenyl)-	
	xanthylium, monochloride	
Synonyms:	RH110, Rhodamine 560, Rhodamine N	HO
MF:	$C_{20}H_{15}N_2O_3 \bullet CI$	
FW:	366.8	
Purity:	≥97%	
Ex./Em. Max:	496/520 nm	
Supplied as:	A crystalline solid	H ₂ N NH ₂
Storage:	-20°C	• CI ⁻
Stability:	≥4 years	
		•

Special Conditions: Keep cool and dry. Protect from light and moisture. Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Rhodamine 110 (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the rhodamine 110 (chloride) in the solvent of choice, which should be purged with an inert gas. Rhodamine 110 (chloride) is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide. Rhodamine 110 (chloride) is slightly soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Rhodamine 110 is a green fluorescent cationic dye with excitation and emission maxima of 496 and 520 nm, respectively.¹ When incorporated with a hydrolytic substrate (e.g., proteinase or peptidase substrates), it can be used as a highly sensitive detection reagent in fluorescence-based enzyme assays.² Rhodamine 110 has also been used in a fluorescence quenching method for determining trace nitrite and as a probe for cytochrome P450 activity.^{3,4}

References

- 1. Thatte, H.S., Rhee, J.-H., Zagarins, S.E., et al. Acidosis-induced apoptosis in human and porcine heart. Ann. Thorac. Surg. 77(4), 1376-1383 (2004).
- 2. 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).
- 3. Zhang, X., Wang, H., Fu, N.-N., et al. A fluorescence quenching method for the determination of nitrite with Rhodamine 110. Spectrochimica Acta. A. Mol. Biomol. Spectrosc. 59(8), 1667-1772 (2003).
- Yatzeck, M.M., Lavis, L.D., Chao, T.-Y., et al. A highly sensitive fluorogenic probe for cytochrome P450 4. activity in live cells. Bioorg. Med. Chem. Lett. 18(22), 5864-5866 (2008).

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

SAFFTY 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

uyer 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, 10/06/2022

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