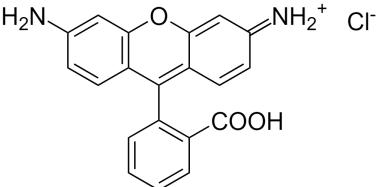


Rhodamine 110 *CAS 13558-31-1*Catalog number: 86
Unit size: 1 g**Product Details**

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

Chemical Properties

Appearance	Red solid solid
Molecular Weight	366.8
Soluble In	DMSO
Chemical Structure	 The chemical structure of Rhodamine 110 is shown. It features a central carbon atom bonded to a 4-aminophenyl group, a 4-carboxyphenyl group, and two 4-aminophenyl groups. One of the 4-aminophenyl groups is further substituted with a chloride ion (Cl-). The structure is a derivative of a diarylmethane, where the two phenyl rings are substituted with amino groups and a central methine group is substituted with a carboxyphenyl group.

Spectral Properties

Excitation Wavelength	500 nm
Emission Wavelength	522 nm

Applications

Rhodamine 110 (R110) is the precursor for a variety of R110-based enzyme substrates. It is also used as a calibration standard for R110-based enzyme substrates. Compared to AMC and AFC enzyme substrates, fluorogenic R110 substrates are generally more sensitive, giving longer absorption and fluorescence wavelength than AMC and AFC substrates.