

Ubiquitin-Rhodamine 110 Substrate

Catalog Number: 12901

Unit Size: 50 ug

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--

Expiration Date	24 months upon receiving
-----------------	--------------------------

Chemical Properties

Appearance	Solid
------------	-------

Molecular Weight	8933.9
------------------	--------

Soluble In	DMSO
------------	------

Spectral Properties

Excitation Wavelength	500 nm
-----------------------	--------

Emission Wavelength	522 nm
---------------------	--------

Applications

Ubiquitin-Rhodamine 110 Substrate consists of full-length ubiquitin conjugated to Rhodamine 110. In its intact form, the substrate exhibits minimal fluorescence. Cleavage by deubiquitinating enzymes releases fluorescent Rhodamine 110, producing a measurable green signal with excitation and emission maxima of approximately 500 nm and 522 nm. This format enables real-time and quantitative measurement of DUB activity.

The substrate is commonly used in biochemical assays for characterizing DUB activity, determining kinetic parameters, and screening potential modulators. Its fluorescence properties are compatible with standard fluorescence microplate readers and high-throughput assay platforms.