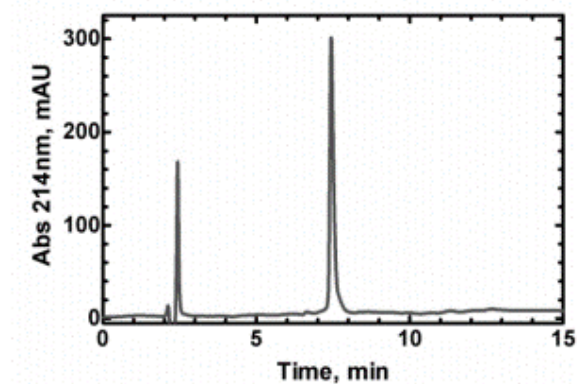


Product Information

Description:	Ubiquitin-rhodamine 110 is a quenched, fluorescent substrate for deubiquitylases, especially ubiquitin C-terminal hydrolases. Cleavage of the amide bond between the C-terminal glycine of ubiquitin and rhodamine results in an increase in rhodamine fluorescence at 535 nm (Exc. 485 nm).
Supplied as:	Lyophilized Solid
Solubility:	Aqueous buffers, DMSO
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	8934.2 Da
Purity:	≥95% determined by RP-HPLC
Applications:	Measurement of deubiquitylase activity.

Quality Control Data

RP-HPLC



Deconvoluted mass spectrum

