

## Ubiquitin- $\epsilon$ -Lys-TAMRA-Gly-OH , DUB Fluorescence Polarization Substrate

Cat. # SI270

**Background:**

Fluorescence polarization substrate for deubiquitinating enzymes. The substrate consists of a 5-carboxytetramethylrhodamine (TAMRA)-labeled Lys-Gly dipeptide where the  $\epsilon$ -amino group of Lys is linked to the C-terminus of ubiquitin. The substrate allows continuous fluorescence polarization measurement of cleavage of the isopeptide bond between ubiquitin and the  $\epsilon$ -amino group of Lys. The rhodamine fluorophore minimizes interference from compounds and assay components that fluoresce at lower wavelengths.

**Application:**

- Determining the activity and specificity of isopeptidases for Ub/Ubl-proteins
- High Throughput Screening for modulators of DUB activity

**Product Information**

<b>Purity:</b>	$\geq$ 95% by RP-HPLC
<b>Molecular Weight:</b>	9162 Da by MS (calc MW: 9165 Da)
<b>Physical State:</b>	Dry powder
<b>Quantity:</b>	50 $\mu$ g
<b>Solubility:</b>	DMSO.
<b>Storage:</b>	-80°C. Avoid repeated freeze/thaw cycles. Protect from light.
<b>Typical assay concentration:</b>	10-100nM

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2009 LifeSensors, Inc. All Rights Reserved