

USP8

(Ubiquitin carboxyl-terminal hydrolase 8, Deubiquitinating enzyme 8, Ubiquitin isopeptidase Y, Ubiquitin thiolesterase 8, Ubiquitin-specific-processing protease 8, KIAA0055, UBPY, UBP8)

- **CAT#:** USP8
- **Enzyme:** Human USP8
- **Substrate:** Fluorogenic substrate [Ubiquitin-AMC], 100 nM
- **Reference Compound IC50s:**

Compounds	IC50 (μ M)
UCHL1 I	8.3
UIPase I	>200
UI1	72.9

- **Scientific Information:** P40818 (<http://www.expasy.org/uniprot/P40818>)

Reference Compound IC50 for USP8

