

USP14

(Ubiquitin carboxyl-terminal hydrolase 14, Deubiquitinating enzyme 14, Ubiquitin thiolesterase 14, Ubiquitin-specific-processing protease 14, TGT, UBP14)

- **CAT#:** USP14
- **Enzyme:** Human USP14
- **Substrate:** Fluorogenic substrate [Ubiquitin-AMC], 100 nM
- **Reference Compound IC50s:**

Compounds	IC50 (μ M)
UCHL1 I	72.4
Ub Aldehyde	2.8
UCHL1 2	76.5

- **Scientific Information:** P54578 (<http://www.expasy.org/uniprot/P54578>)

Reference Compound IC50 for USP14

