

BAP1

(Ubiquitin carboxyl-terminal hydrolase BAP1, BRCA1-associated protein 1, Cerebral protein 6, KIAA0272, hucep-6)

- **CAT#:** BAP1
- **Enzyme:** Human BAP1
- **Substrate:** Fluorogenic substrate [Ubiquitin-AMC], 100 nM
- **Reference Compound IC50s:**

Compounds	IC50 (μ M)
UCHL1 I	180
UIPase I	>300
UI1	>300

- **Scientific Information:** Q92560 (<http://www.expasy.org/uniprot/Q92560>)

Reference Compound IC50 for BAP1

