

## Product Information

---

<b>Description:</b>	BAP1 is a deubiquitinase that plays a key role in chromatin by mediating deubiquitination of histone H2A and HCFC1. It acts as a regulator of cell growth.
<b>Background:</b>	BAP1 plays a critical role in regulating various cellular processes, including DNA repair, cell cycle control, and gene expression. It is particularly known for its involvement in maintaining genomic stability by deubiquitinating histones and other proteins involved in chromatin remodeling. Mutations or loss of BAP1 function have been associated with a higher risk of developing certain cancers, including uveal melanoma and mesothelioma, underscoring its significance as a guardian of genome integrity.
<b>Tag:</b>	His-Tag
<b>Species:</b>	Human
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥85% by SDS-PAGE
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>MW:</b>	82 kDa
<b>Activity:</b>	This enzyme cleaves Ub-AMC.