

Product Information

Description:	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.
Background:	BAP1 is the first large, nuclear-localized UCH to be identified. It is a protein of 90 kDa that binds to the BRCA RING finger and enhances BRCA1-mediated cell growth suppression.
Tag:	His-Tag
Species:	Human
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	82 kDa
Activity:	This enzyme cleaves Ub-AMC.