

## CUTANA™ Protease Inhibitor Tablets

<b>Catalog No</b>	21-1027-S	<b>Pack Size</b>	25 Small Tablets
<b>Lot No</b>	24292001-81	<b>25X Stock</b>	420 µL per tablet

### DESCRIPTION

This EDTA-free cocktail inhibits various serine and cysteine proteases, resulting in stabilization of target proteins isolated from animal tissues, plants, and yeast for *in situ* assays like ChIC/CUT&RUN (EpiCypher 14-1048) or CUT&Tag (EpiCypher 14-1102/14-1103). Absence of EDTA and other chelating agents is essential for the controlled activation of pAG-MNase (EpiCypher 15-1016/15-1116) and pAG-Tn5 (EpiCypher 15-1017/15-1117), but leaves metallo- and aspartic proteases inhibited. This product is also used in the preparation of CUTANA™ Nuclei Extraction Buffer (EpiCypher 21-1026).

### TECHNICAL INFORMATION

<b>Storage</b>	Tablets stored in their original container are stable at 4°C until the date of expiration on the label.		
<b>Formulation</b>	Proprietary mixture of protease inhibitors; EDTA free.		
<b>Instructions for Use</b>	Reconstitute each tablet in 420 µL molecular biology-grade water to prepare a 25X stock. Leftover 25X protease inhibitor can be stored for 12 weeks at -20°C.		
<b>Reactions/Tablet</b>	<b>CUT&amp;RUN</b>	~5 reactions	
	<b>CUT&amp;Tag</b>	~6 reactions	
	<b>Nuclei Extraction</b>	~40 reactions	