

CUTANA™ Protease Inhibitor Tablets

Catalog No	21-1027-L	Pack Size	20 Large Tablets
Lot No	24292001-82	25X Stock	2 mL 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

Storage	Tablets stored in their original container are stable at 4°C until the date of expiration on the label.		
Formulation	Proprietary mixture of protease inhibitors; EDTA free.		
Instructions for Use	Reconstitute each tablet in 2 mL molecular biology-grade water to prepare a 25X stock. Leftover 25X protease inhibitor can be stored for 12 weeks at -20°C.		
Reactions/Tablet	CUT&RUN	~27 reactions	
	CUT&Tag	~30 reactions	
	Nuclei Extraction	~204 reactions	