

## CUTANA™ Protease Inhibitor Tablets

|                   |             |                  |                   |
|-------------------|-------------|------------------|-------------------|
| <b>Catalog No</b> | 21-1027-S   | <b>Pack Size</b> | 25 Small Tablets  |
| <b>Lot No</b>     | 25038002-81 | <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 |  |