

## CUTANA™ Uncharged pAG-Tn5 for CUT&Tag

<b>Catalog No</b>	15-1025	<b>Species</b>	<i>E. coli</i>
<b>Lot No</b>	20142001-R1	<b>Source</b>	<i>E. coli</i>
<b>Pack Size</b>	20 µL	<b>Epitope Tag</b>	None
<b>Concentration</b>	5.5 µM (dimer)	<b>MW</b>	157 kDa (dimer)

### DESCRIPTION

Products in EpiCypher's IDEA Toolbox (Innovation and Discovery of Epigenetic Applications) offer access to reagents without known or fully defined uses, enabling researchers to explore cutting-edge applications. Due to their novelty and unexplored potential, EpiCypher will engage in limited technical support.

CUTANA Uncharged pAG-Tn5 is a fusion of Proteins A and G to *E. coli* transposase (Tn5) and is the key enzyme for CUT&Tag [1]. This product is highly purified to remove contaminating *E. coli* DNA. For superior normalization, as well as antibody validation and reaction monitoring, SNAP-CUTANA™ nucleosome spike-ins (e.g. EpiCypher 19-1002) are recommended. **This Tn5 is uncharged and must be loaded with user-designed mosaic adapter DNA prior to use in CUT&Tag.** The uncharged protein exists as a monomer but will dimerize when charged with DNA. For a charged enzyme that is ready to use, see EpiCypher 15-1017.

### RECOMMENDED ACCESSORY REAGENTS

<u>Item</u>	<u>CAT</u>	<u>Item</u>	<u>CAT</u>
Anti-Rabbit Secondary Antibody	13-0047	H3K4me3 Positive Ctrl Antibody	13-0041
Anti-Mouse Secondary Antibody	13-0048	Rabbit IgG Negative Ctrl Antibody	13-0042
Magnetic Separation Rack, 0.2 mL	10-0008	CUTANA™ ConA Beads	21-1401
Magnetic Separation Rack, 1.5 mL	10-0012	SNAP-CUTANA™ K-MetStat	19-1002
CUT&RUN 8-strip Tubes	10-0009	Non-HS 2X PCR Master Mix	15-1018

### TECHNICAL INFORMATION

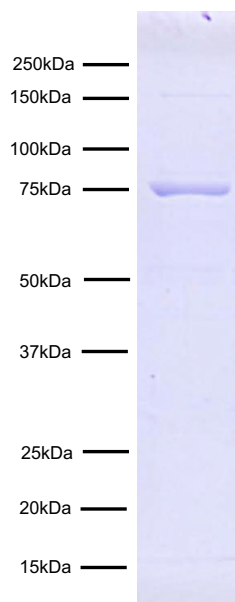
<b>Storage</b>	Stable for one year at -20°C from date of receipt. The protein is not subject to freeze/thaw under these conditions.
<b>Formulation</b>	50 mM HEPES-KOH pH 7.2, 100 mM NaCl, 0.1 mM EDTA, 1 mM DTT, 0.1% Triton X-100, 50% glycerol

### APPLICATION NOTES

pAG-Tn5 transposomes can be assembled as previously described [1]. In brief, 3.2 µL of an equimolar mixture of pre-annealed user-defined Adapter-A and user-defined Adapter-B oligonucleotides (50 µM each, 100 µM total adapter DNA) should be mixed with 20 µL of 5.5 µM pAG-Tn5 fusion protein dimer (a 3:1 molar ratio of adapter DNA to pAG-Tn5 dimer). The mixture is then incubated on a rotating platform for 1 hour at room temperature and stored at -20°C. Specific activity definition of the charged pAG-Tn5 is highly recommended before use in CUT&Tag. **Due to the confounding variable of user-supplied mosaic adapters, EpiCypher will not engage in protocol troubleshooting for this reagent.** For a pre-charged enzyme that is eligible for technical support, see EpiCypher 15-1017.

### REFERENCES

[1] Kaya-Okur et al. *Nat. Commun.* (2019). PMID: 31036827



**FIGURE 1 Protein gel data.** CUTANA pAG-Tn5, Uncharged (1  $\mu$ g) was resolved via SDS-PAGE and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.

US Pat. No. 10,689,643, 11,306,307, EU Pat. No. 3,688,157, 2,999,784  
and related patents and pending applications