

## pA-Tn5 transposase (loaded)

Cat. No. **C01070001**

Format: 15 µl (32 rxns) / 30 µl (64 rxns)

Molecular weight: 67 kD

### Product description

**pA-Tn5 Transposase** is a fusion protein of hyperactive Tn5 transposase and protein A developed for the **CUT&Tag** assay. For ease of use, the fusion protein is pre-loaded with sequencing adapters suitable for single or dual indexing in single or paired-end Illumina platforms. The adaptors contain 19-mer Tn5 mosaic ends and the sequences for PCR amplification with barcoded i7/i5.

Mosaic end\_reverse: [PHO]CTGTCTCTTATACACATCT

Mosaic end\_Adapter A: **TCGTCGGCAGCGTC**AGATGTGTATAAGAGACAG

Mosaic end\_Adapter B: **GTCTCGTGGGCTCGG**AGATGTGTATAAGAGACAG

Underlined regions correspond to the double-stranded part of the adapter, recognized by the transposase. **Bold regions** represent sequences required per PCR amplification with barcoded i7/i5 primers as described by Buenrostro et al (2015).

The protein A has a high affinity to rabbit polyclonal antibodies, mouse IgG2a, IgG2b and IgA, guinea pig IgG, dog IgG, pig IgG. However, the use of secondary antibody (e.g. guinea pig anti-rabbit) is recommended for a higher sensitivity of **CUT&Tag** assay.

### Suggested dilution:

For the standard **CUT&Tag assay**, the recommended dilution of **pA-Tn5 Transposase (loaded)** is 1:250 (0.4 µl of pA-Tn5 for 100 µl of buffer). Please note that depending on the starting amount of cells and/or primary antibody, different dilution in a range 1:50-1:500 might be tested.

### Storage conditions

Store at -20°C.

### Storage buffer

Supplied in solution containing 50% v/v glycerol.

### Precautions

This product is for research use only. Not for use in diagnostic or therapeutic procedures.

#### Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK  
Rue du Bois Saint-Jean, 3  
4102 Seraing - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

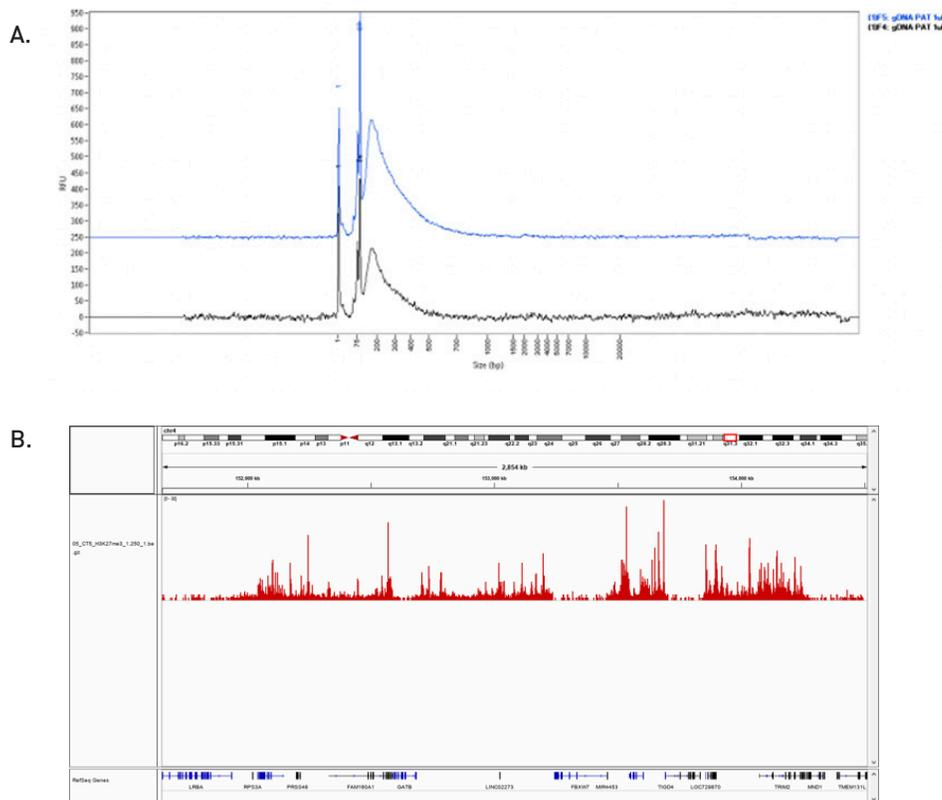
#### Diagenode LLC. USA | NORTH AMERICA

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

Last update: August, 2022

## Quality control

Each lot of pA-Tn5 transposase is quality checked by an in vitro activity test (cleavage of human genomic DNA) (Figure 1, A) and by a CUT&Tag assay using H3K27me3 polyclonal ChIP-seq grade antibody (Cat. No. C15410195) (Figure 1, B).



**Figure 1: Quality control of pA-Tn5 transposase loaded with sequencing adapters**

A: The Fragment Analyzer trace showing the representative cleavage pattern of gDNA. The pA-Tn5 fusion protein (Cat. No. C01070001) efficiently digests gDNA to a smear. 500 ng of human genomic DNA were incubated for 7 min at 55°C with 1 µl of pA-Tn5 fusion protein loaded with appropriated adaptors in a tagmentation buffer (40mM Tris-HCl pH7.5, 40mM MgCl<sub>2</sub> and 12.5% DMF). The reaction was stopped by adding SDS, cleaned-up and resolved on the Fragment Analyzer to assess the cleavage.

B: Representative screenshot at selected locus obtained using Diagenode pA-Tn5 fusion protein (Cat. No. C01070001) and H3K27me3 polyclonal ChIP-seq grade antibody (Cat. No. C15410195) following CUT&Tag protocol (Kaya-Okur, H.S., Nat Commun 10, 1930 [2019]).

### Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK  
Rue du Bois Saint-Jean, 3  
4102 Seraing - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

### Diagenode LLC. USA | NORTH AMERICA

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com