

Tagmentation Test Protocol for Tn5 or pA-Tn5

You will need:

- Assembled Tn5 or pA-Tn5 Transposomes
- Tagmentation Reaction Buffer (appropriate for your assay)
- Genomic DNA, 50 ng/ μ L
- Freshly prepared 0.5% SDS
- Thermal cycler
- Columns or beads for purification
- Agilent TapeStation or similar

Prepare your Assembled Tn5 or pA-Tn5 Transposomes, and Tagmentation Reaction Buffer appropriate for your assay.

1. Prepare a mastermix with your Tagmentation Reaction Buffer and genomic DNA (keep genomic DNA constant at 150 ng, and titrate volume of Assembled Transposomes).
2. Aliquot each sample and add remaining water to bring to volume (25 μ L minus the Assembled Transposomes volume).
3. Add Assembled Transposomes (for example, titrate 1 μ L, 2 μ L, 4 μ L).
4. Mix the reaction and leave at 55°C for 7 minutes.
5. Add 2.5 μ L 0.5% SDS and incubate for another 3 minutes at 55°C.
6. Column or bead purify and elute in 10 μ L 10 mM TE.
7. Run on Agilent TapeStation.

Technical Support

If you need assistance at any time, please call or send an e-mail to Active Motif Technical Service at one of the locations listed below.

Active Motif North America

1914 Palomar Oaks Way, Suite 150
Carlsbad, CA 92008, USA

Toll Free: 877 222 9543
Direct: 760 431 1263
Fax: 760 431 1351
E-mail: tech_service@activemotif.com

Active Motif Europe

Waterloo Atrium
Drève Richelle 167 – boîte 4
BE-1410 Waterloo, Belgium

Germany Free Phone: 0800 181 99 10
France Free Phone: 0800 90 99 79
UK Free Phone: 0800 169 31 47
Fax: +32 (0)2 653 0050
E-mail: eurotech@activemotif.com

Active Motif Japan

Azuma Bldg, 7th Floor
2-21 Ageba-Cho, Shinjuku-Ku
Tokyo, 162-0824, Japan

Direct: +81 3 5225 3638
Fax: +81 3 5261 8733
E-mail: japantech@activemotif.com

Active Motif China

889 Wanhangu Road
Suite 1602B
Jingan District
Shanghai, 200040, China

Direct: (86)-21-20926090
Cell Phone: 18521362870
E-mail: techchina@activemotif.com

Visit Active Motif online at <http://www.activemotif.com>

For Research Use Only. Not for use in diagnostic procedures.

