

pAM_1C_JunD Vector

Catalog No.: 53044

Format: 50 µg

Description

The pAM_1C_JunD Vector is designed for use with Active Motif's Tag-ChIP-IT® Kit (Catalog No. 53022). The DNA sequence for transcription factor JunD (NCBI Accession NM_005354.5) was cloned into the pAM_1C empty vector using InFusion Cloning® (Clontech) to append the AM-tag sequence to the carboxy-terminus. pAM_1C_JunD contains the Human Beta-actin promoter (ACTB) which provides constitutive high-level protein expression in mammalian systems. The vector may be used for transient transfections, or puromycin selection can be applied to select for stable gene expression. Western blot analysis with the AM-Tag antibody (Catalog No. 61677) can confirm tagged-protein expression. Cells expressing the fusion protein are ready for use in the Tag-ChIP-IT Kit.

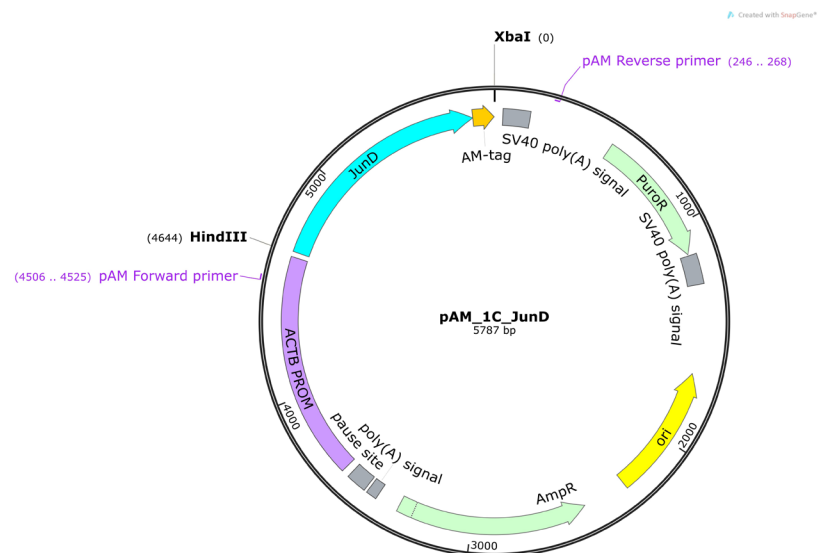
Contents

- 50 µg of pAM_1C_JunD Vector provided as lyophilized DNA.
Resuspend plasmid to a concentration of 1 µg/µl in sterile H₂O. Store reconstituted DNA at -20°C.

pAM_1C_JunD Vector Features and Circle Map

The following features are present in the pAM_1C_JunD Vector based on nucleotide sequence.

SV40 poly(A) signal	39-160
pAM Reverse primer	246-268
Puromycin coding region	542-1141
SV40 poly(A) signal	1148-1282
Col/E1-derived plasmid origin of replication	1685-2273
β-lactamase (Amp ^r) coding region	2473-3333
Synthetic polyadenylation signal	3438-3486
RNA pol II transcriptional pause site	3500-3591
ACTB promoter	3626-4626
pAM Forward primer	4505-4525
JunD	4650-5690
AM-tag	5691-5786



Quality Control

Plasmid construct has been confirmed by restriction analysis and sequence verified.

Shipping & Storage

Products are shipped at room temperature.

Lyophilized DNA is stable for 12 months when stored at -20°C.

Resuspended DNA is stable for 6 months when stored at -20°C. Avoid repeated freeze/thaw cycles.

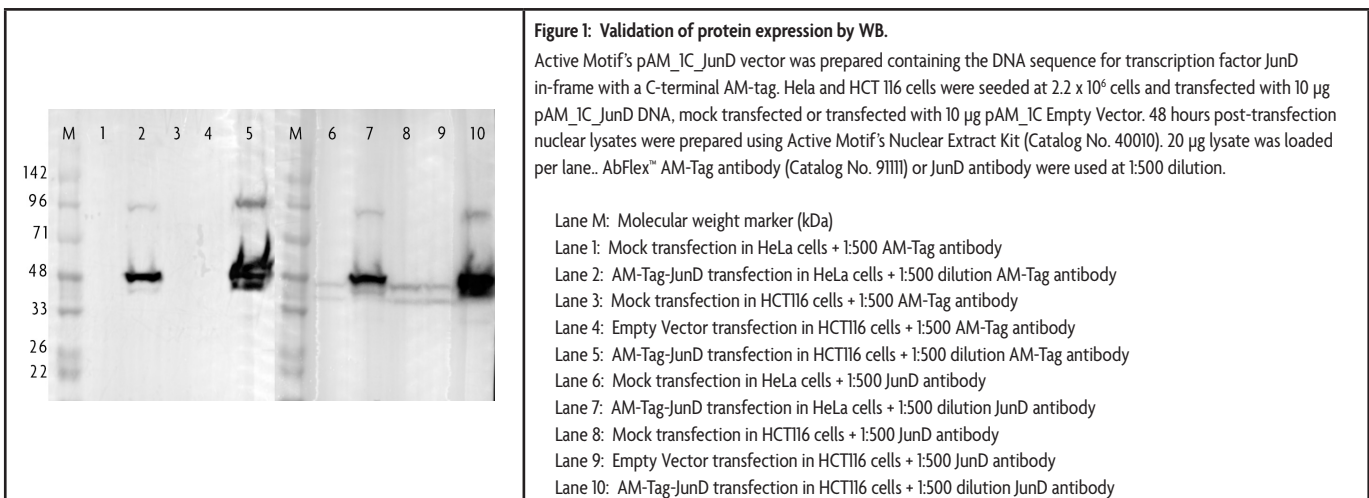
Background

JunD is a component of the activator protein-1 (AP-1) transcription factor family. AP-1 proteins play a role in the expression of many genes involved in proliferation and cell cycle progression including neuronal apoptosis, learning process, drug-induced behavioral responses, bone growth and differentiation, and embryo development. AP-1 is composed of a mixture of heterodimeric complexes of proteins derived from the Fos and Jun families including c-Fos, FosB, Fra-1, Fra-2, c-Jun, JunB and JunD. Only Jun proteins can form transcriptionally active homodimers. The N-terminus of Jun proteins is involved in transcriptional activation, while the C-terminus is involved in dimerization and DNA binding. AP-1 dimers bind the TPA-response element (TRE). JunD exists as two distinct isoforms (40 kDa & 45 kDa) generated by the use of two translational start sites within the JunD mRNA. Full length JunD contains an additional 48 amino acids at the amino-terminus. Both isoforms are expressed in all cell types at approximately the same levels.

GENERAL PRODUCT USE

pAM_1C_JunD can be used as a control vector in Active Motif's Tag-ChIP-IT® Kit (Catalog No. 53022). We have validated this vector using the following transfection conditions in HCT116 cells (human colorectal carcinoma). Optimization of transfection conditions may be required for use in other cell types.

1. In a 100 mm dish, seed 2.2×10^6 HCT116 cells per dish in McCoy's 5A medium supplemented with 10% fetal bovine serum. Incubate in a humidified incubator for 24 hours. We recommend setting up two dishes. One dish can be used for chromatin preparation and the second dish can be used to confirm expression of the tagged protein by Western blot.
2. Prepare a separate microcentrifuge tube for each transfection reaction. Add 10 μ g pAM_1C_JunD construct to Opti-MEM media in a final volume of 550 μ l.
3. Add 30 μ l FuGENE HD Transfection Reagent (Catalog No. 32042) drop wise directly to the media/DNA mixture. Do not allow FuGENE to come directly into contact with the plastic. Mix the solution by pipetting up and down and incubate at room temperature for 15-30 minutes.
4. Add 580 μ l of the media/DNA/FuGENE mixture drop wise to each 100 mm dish. Gently swirl dish or incubate on a shaker at 100 rpm for 2 minutes to evenly distribute transfection mixture.
5. Return dish to humidified incubator for 48 hours before proceeding with the chromatin preparation.
6. Refer to the Tag-ChIP-IT manual to prepare chromatin.
7. Use Active Motif's Nuclear Extract Kit (Catalog No. 40010) to prepare nuclear lysates from the second dish for detection of the tagged protein as compared to untransfected cells. Use 20 μ g nuclear lysate denatured at 98C for 10 minutes in sample loading buffer per well of a 4-20% Tris-Glycine gel. Run at 90 mAmps until the dye front runs off the gel. Transfer to a nitrocellulose membrane for 90 minutes at 35 volts. Block the membrane with 5% non-fat milk in PBS for 30 minutes. Add the AbFlex™ AM-Tag antibody (Catalog No. 91111) at a 1:500 - 1:1000 dilution in 5% non-fat milk in PBS for 1 hour. Wash the membrane three times for 5 minutes with distilled water. Add an anti-mouse HRP-conjugated secondary at 1:1000 dilution for 1 hour in 5% non-fat milk in PBS. Wash the membrane three times for 5 minutes with PBS containing 0.05% Tween 20 followed by 3-5 rinses of the membrane with distilled water. Detect with ECL or chemiluminescent substrate.



Technical Services

If you need assistance at any time, please call Active Motif Technical Service at one of the numbers listed below.

Active Motif North America

1914 Palomar Oaks Way, Suite 150
Carlsbad, CA 92008

USA

Toll Free: 877 222 9543

Telephone: 760 431 1263

Fax: 760 431 1351

E-mail: tech_service@activemotif.com

Active Motif Europe

Avenue Reine Astrid, 92

B-1330 La Hulpe, Belgium

UK Free Phone: 0800 169 31 47

France Free Phone: 0800 90 99 79

Germany Free Phone: 0800 181 99 10

Telephone: +32 (0)2 653 0001

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

Telephone: +81 3 5225 3638

Fax: +81 3 5261 8733

E-mail: japantech@activemotif.com

Active Motif China

787 Kangqiao Road

Building 10, Suite 202, Pudong District

Shanghai, 201315, China

Telephone: (86)-21-20926090

Hotline: 400-018-8123

Visit Active Motif on the worldwide web at <http://www.activemotif.com>

At this site:

- Read about who we are, where we are, and what we do
- Review data supporting our products and the latest updates
- Enter your name into our mailing list to receive our catalog, *MotifVations* newsletter and notification of our upcoming products
- Share your ideas and results with us
- View our job opportunities

Don't forget to bookmark our site for easy reference!