

# Chariot™

Simple, efficient protein delivery

(version C)

Catalog Nos. 30025 & 30100

## Active Motif North America

1914 Palomar Oaks Way, Suite 150

Carlsbad, California 92008, USA

Toll free: 877 222 9543

Telephone: 760 431 1263

Fax: 760 431 1351

## Active Motif Europe

104 Avenue Franklin Roosevelt

B-1330 Rixensart, 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

## 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

TABLE OF CONTENTS	Page
<b>Overview</b>	
Description .....	1
Kit Contents .....	1
Quality Control .....	1
<b>Introduction</b> .....	2
<b>Flow Chart of Process</b> .....	3
<b>Protocols</b>	
Preparation of Chariot .....	4
Preparation of $\beta$ -galactosidase .....	4
Transfection Protocol .....	4
Transfection Protocol for Suspension Cells .....	7
$\beta$ -galactosidase Staining Protocol .....	8
<b>References</b> .....	9
<b>Troubleshooting Guide</b> .....	10
<b>Related Products</b> .....	11
<b>Technical Services</b> .....	12

## Overview

Chariot™ is a patented\* transfection reagent suitable for the delivery of proteins, peptides and antibodies into a variety of cultured mammalian cells. Chariot Kits come in two sizes and can be used with many different plate sizes as indicated in the table below.

Plate size	Chariot (μl)	No. of Transfections per Kit (25-reaction Kit)	No. of Transfections per Kit (100-reaction Kit)
96-well	0.5	300	1200
24-well	2	75	300
12-well	3.5	42	171
6-well or 35 mm	6	25	100
60 mm	20	7	30
100 mm	50	3	12

## Kit Contents

Component	Quantity		Composition	Storage
	25 rxns	100 rxns		
Chariot Transfection Reagent	0.3 mg	1.2 mg		-20°C for up to 6 months
PBS	1 ml	1 ml	1.5 mM KH <sub>2</sub> PO <sub>4</sub> 150 mM NaCl 5 mM Na <sub>2</sub> HPO <sub>4</sub>	-20°C
β-galactosidase (Positive Control)	25 μg	25 μg		-20°C for up to 6 months
Sterile H <sub>2</sub> O	1 ml	1 ml		-20°C

## Quality Control

- Chariot Transfection Reagent is functionally tested by delivery of protein and antibody into many different cell lines (including HS-68, NIH 3T3 and HeLa). Please visit our website for a complete list of cell lines that have been tested.
- Chariot Transfection Reagent is tested for the absence of bacterial and fungal contamination in DMEM cell culture media supplemented with 10% FBS.
- Chariot Transfection Reagent is non-cytotoxic at the recommended concentrations.

\* Chariot is covered under U.S. Patent No. 6,841,535. Purchase includes the right to use for basic research purposes only. Other use licenses are available, please contact Active Motif Technical Services for additional information.

## Introduction

---

Chariot™ is a revolutionary new transfection reagent capable of efficiently delivering proteins, peptides and antibodies into cultured mammalian cells in less than two hours.

Current transfection techniques include microinjection<sup>1</sup>, calcium phosphate coprecipitation<sup>2</sup>, cationic liposomes<sup>3</sup>, viral vectors<sup>4</sup> and electroporation<sup>5</sup>. These methods are capable of transporting DNA into cells, but the techniques can be cumbersome and even cytotoxic. Once transfection has been completed, the researcher must wait 12-80 hours post-transfection to detect expression of the gene of interest.

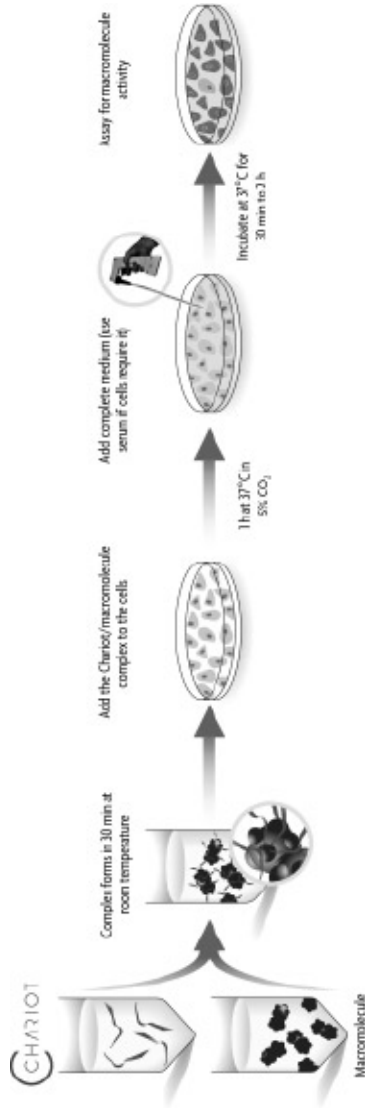
Recently, novel methods to deliver proteins have been reported and commercialized, *i.e.* HIV-1 TAT, *Drosophila* Antennapedia homeotic transcription factor and herpes simplex virus-1 DNA binding protein VP22 (for a recent review, see ref. 6). Transduction via TAT-fusion proteins results in inactivation and denaturation of the protein. To deliver an active protein, correct renaturation is required upon internalization. In addition, TAT must be covalently linked to the compound or macromolecule to be delivered by a chemical reaction.

Penetratin 1 is a 16 amino acid peptide corresponding to the third helix of the homeodomain of Antennapedia protein. Activated Penetratin has an N-terminal pyridyl disulfide that is used to covalently couple it to the macromolecule to be delivered. However, chemical coupling can be cumbersome and requires the macromolecule being delivered to carry a free thiol group.

Another delivery system utilizes the translocation properties of the 38 kDa herpes simplex virus-1 DNA binding protein VP22. VP22 must be fused to the peptide/protein to be delivered and, therefore, requires the construction of a suitable expression vector.

Chariot is a 2843 dalton peptide and unlike these current transfection techniques, Chariot forms a non-covalent complex with the protein, peptide or antibody of interest. This completely bypasses the transcription-translation process associated with gene expression, reducing the time until the cells can be assayed from days to under two hours. The Chariot-macromolecule complex stabilizes the macromolecule and helps to protect it from degradation during the transfection process<sup>7, 8</sup>. Upon internalization, the complex dissociates and the macromolecule is free to proceed to its target organelle. Moreover, efficient delivery is observed at 4°C, suggesting that the delivery mechanism is independent of the endosomal pathway<sup>8</sup>. Therefore, the macromolecule is not subjected to the conditions of that pathway, which can modify the structure of the macromolecule during internalization. Chariot is non-cytotoxic and serum-independent<sup>8</sup>. It is ideal for *in vivo* studies because fixing is not required.

# Flow Chart of Process



## Protocols

---

### Preparation of Chariot

Chariot is provided as a lyophilized powder and arrives at room temperature. Store at  $-20^{\circ}\text{C}$  for long term storage.

25-reaction Kit: 0.3 mg of Chariot is provided. Resuspend the entire contents of the Chariot vial in 150  $\mu\text{l}$  sterile  $\text{H}_2\text{O}$ . Mix gently by tapping tube.

100-reaction Kit: 1.2 mg of Chariot is provided. Resuspend the entire contents of the Chariot vial in 600  $\mu\text{l}$  sterile  $\text{H}_2\text{O}$ . Mix gently by tapping tube.

After resuspension, store at  $-20^{\circ}\text{C}$  in small aliquots to avoid repeated freeze/thaw cycles.

**NOTE:** If you will be transfecting a peptide or low molecular weight protein ( $< 10$  kDa), you must make an additional 1:10 dilution of the resuspended Chariot solution in sterile  $\text{H}_2\text{O}$  before use. This will improve transfection efficiency by preventing the formation of aggregates.

**NOTE:** Diluting Chariot in DMSO has been shown to improve delivery of some antibodies. It has also provided more uniform delivery of some proteins. Please refer to the sections in the Troubleshooting Guide (page 10) for more information.

### Preparation of $\beta$ -galactosidase

$\beta$ -galactosidase is an enzyme that hydrolyses  $\beta$ -galactosides, such as lactose and the artificial chromogen X-gal.  $\beta$ -galactosidase is composed of four identical subunits. When  $\beta$ -galactosidase hydrolyses X-gal, it produces a blue color that can be visualized under a bright field microscope. A 119 kDa subunit of the  $\beta$ -galactosidase protein is provided in this kit as a positive control for the Chariot Transfection Reagent.

The  $\beta$ -galactosidase positive control protein is provided as a lyophilized powder. Store at  $-20^{\circ}\text{C}$ . 25  $\mu\text{g}$  of protein is provided. Resuspend the entire contents of the  $\beta$ -galactosidase positive control vial in 100  $\mu\text{l}$  sterile  $\text{H}_2\text{O}$ . This makes a 0.25  $\mu\text{g}/\mu\text{l}$  stock. Use 1.0  $\mu\text{g}$   $\beta$ -galactosidase per transfection reaction. After resuspension, aliquot and store at  $-20^{\circ}\text{C}$  to avoid repeated freeze/thaw cycles.

When the  $\beta$ -galactosidase positive control protein is transfected, the cells must be stained to assay for the efficiency of the transfection. A detailed staining protocol and buffer recipes are provided on page 8. For your convenience, Active Motif also offers a  $\beta$ -galactosidase Staining Kit (Catalog No. 35001) available separately.

## Transfection Protocol

**These conditions are recommended as guidelines only.** Efficient transfection may require optimization of reagent concentration, cell number and exposure time of cells to the Chariot-macromolecule complex. Conditions should also be optimized for each cell line and kept consistent to obtain reproducible results. This procedure has been optimized for the transfection of adherent cells using the  $\beta$ -galactosidase positive control. A protocol for using suspension cells can be found on page 7.

1. Seed cells according to the following table:

Plate	Surface Area (mm <sup>2</sup> )	Cells	Growth Medium (ml)
96-well	20	1.0 x 10 <sup>4</sup>	0.2
24-well	200	0.05 x 10 <sup>6</sup>	0.5-1.0
12-well	401	0.1 x 10 <sup>6</sup>	1-2
6-well or 35 mm	962	0.3 x 10 <sup>6</sup>	3-5
60 mm	2827	0.8 x 10 <sup>6</sup>	5
100 mm	7854	2.2 x 10 <sup>6</sup>	10

2. Incubate the cells at 37°C in a humidified atmosphere containing 5% CO<sub>2</sub> until the cells are 40-50% confluent.

**Note:** The transfection efficiency may be sensitive to culture confluency, so it may be necessary to optimize cell density for each cell line.

3. **Preparation of diluted macromolecule:** Dilute your protein, antibody or peptide of interest in the appropriate volume of PBS as indicated in the table below. For antibodies or proteins, we recommend using 0.5-1  $\mu$ g per transfection reaction. For peptides or proteins that are < 10 kDa, we recommend using 100-500 ng per transfection reaction. **NOTE:** The amount of macromolecule used must be optimized, we recommend performing a titration of the protein/antibody or peptide quantities.

Plate	Final volume of macromolecule dilution in PBS ( $\mu$ l)
96-well (or 8-well chamber slide)	10*
24-well	50
12-well	100
6-well or 35 mm	100
60 mm	200
100 mm	200

\* If you are using 96-well plates, you may wish to reduce the amount of protein or antibody used to 0.2  $\mu$ g and amount of peptide to 50 ng.

- 4 **Preparation of diluted Chariot:** In a separate tube, dilute the appropriate volume of Chariot into sterile water as indicated in the table below. For low molecular weight proteins and peptides, a 1/10 dilution of the resuspended Chariot must be made before use to ensure optimal transfection.

Plate	Sterile H2O for Chariot dilution (µl)	Protein/Antibody*	Peptide/LMW Proteins**
		Volume of Chariot (µl)	Volume of a 1/10 dilution of Chariot (µl)
96-well	10	0.5	0.5
24-well	50	2	2
12-well	100	3.5	3.5
6-well or 35 mm	100	6	6
60 mm	200	20	20
100 mm	200	50	50

- For multiple transfections, do not use a master mix that exceeds the volume required for 4 transfections (6-well plate size) per tube, as this may cause aggregation. (**NOTE:** This does not apply to peptide transfections)
- \* Detection of fluorescent antibodies in cells requires a large amount of antibody to be delivered. It may be necessary to increase the amount of Chariot and antibody used.
- \*\* Chariot interacts via hydrophobic interactions. Each peptide or protein will have a different hydrophobicity, which is more apparent with lower molecular weight molecules. The amount of Chariot may need to be titrated to determine the optimal amount.

5. **Preparation of Chariot-macromolecule complex:** Add the diluted macromolecule from Step 3 to the diluted Chariot from Step 4 and incubate at room temperature for 30 minutes to allow the Chariot-macromolecule complex to form.
6. **Transfection of cells:** Aspirate medium from cells to be transfected and wash cells once with PBS.
7. Overlay the cells with the Chariot-macromolecule complex. Next, add the appropriate amount of serum-free medium to the plate as indicated in the table below. Gently rock the plate to ensure even delivery.

Plate	Volume of Chariot-Macromolecule Complex (µl)	Serum-Free Medium (µl)	Final Transfection Volume (µl)
96-well	20	80	100
24-well	100	100	200
12-well	200	150	350
6-well or 35 mm	200	400	600
60 mm	400	1600	2000
100 mm	400	4600	5000

8. Incubate at 37°C in a humidified atmosphere containing 5% CO<sub>2</sub> for one hour.
9. Continue to incubate at 37°C in a humidified atmosphere containing 5% CO<sub>2</sub> for 30 minutes to 2 hours. **OPTIONAL:** If your cells are sensitive to serum starvation you can choose to add complete growth medium to the cells. **Do not** remove the Chariot-macromolecule complex. We suggest adding the minimum amount of complete growth medium needed as the presence of serum can interfere with Chariot transfection efficiency.

**Peptide:** 0.5-1 hour

**Protein:** 1-2 hours

**Antibody:** 2 hours

10. Process the cells for observation or detection assays. Cells may be fixed or observed directly.

### Transfection Protocol for Suspension Cells

**These conditions are recommended as guidelines only.** Efficient transfection may require optimization of reagent concentration, cell number and exposure time of cells to the Chariot-macromolecule complex. Conditions should also be optimized for each cell line and kept consistent to obtain reproducible results.

1. Seed cells according to the following table:

Plate	Surface Area (mm <sup>2</sup> )	Cells	Growth Medium (ml)
96-well	20	1.0 x 10 <sup>4</sup>	0.2
24-well	200	0.05 x 10 <sup>6</sup>	0.5-1.0
12-well	401	0.1 x 10 <sup>6</sup>	1-2
6-well or 35 mm	962	0.3 x 10 <sup>6</sup>	3-5
60 mm	2827	0.8 x 10 <sup>6</sup>	5
100 mm	7854	2.2 x 10 <sup>6</sup>	10

2. The Chariot-macromolecule complex is assembled in the same manner as described for adherent cells. See Steps 3-5 on pages 5-6.
3. Collect the suspension cells by centrifugation at 200-400 x g for 5 minutes. Remove the supernatant.
4. Wash the cells twice with 1X PBS.
5. Centrifuge at 200-400 x g for 5 minutes to pellet the cells. Remove the supernatant.
6. Resuspend the cell pellet in the Chariot-macromolecule complex. Add serum-free medium to achieve the Final Transfection Volume. (See Step 7 page 6)
7. Incubate at 37°C in a humidified atmosphere containing 5% CO<sub>2</sub> for one hour.
8. Continue to incubate at 37°C in a humidified atmosphere containing 5% CO<sub>2</sub> for 30 minutes to 2 hours. **OPTIONAL:** If your cells are sensitive to serum starvation you can choose to

add complete growth medium to the cells. **Do not** remove the Chariot-macromolecule complex. We suggest adding the minimum amount of complete growth medium needed as the presence of serum can interfere with Chariot transfection efficiency.

9. Process the cells for observation or detection assays. Cells may be fixed or observed directly.

### $\beta$ -galactosidase Staining Protocol

#### Make stock solutions:

400 mM potassium ferricyanide	Store at -20°C
400 mM potassium ferrocyanide	Store at -20°C
200 mM magnesium chloride	Store at -20°C
20 mg/ml X-gal 5-bromo-4-chloro-3-indolyl- $\beta$ -D-galactopyranoside resuspended in DMF (N, N-Dimethyl-formamide)	Store at -20°C in the dark
10X PBS	Store at room temperature
10X Fixing Solution 20% formaldehyde 2% glutaraldehyde in 10X PBS	Store at -20°C

#### Staining Protocol

**Note:** Use volumes appropriate for the size of the plate or dish. Ensure that the Fixing and Staining Solutions cover the cells.

1. Remove the growth medium from the transfected cells and rinse 3 times with 1X PBS.
2. Add 1X Fixing Solution. (Dilute the 10X stock in sterile water to make a 1X solution.) Incubate at room temperature for 5-10 minutes.
3. Prepare the Staining Solution. **Staining Solution should be made fresh each time.**

250 $\mu$ l 400 mM potassium ferricyanide	(4 mM final)
250 $\mu$ l 400 mM potassium ferrocyanide	(4 mM final)
250 $\mu$ l 200 mM magnesium chloride	(2 mM final)
1.25 ml X-gal (20 mg/ml in DMF)	(1 mg/ml final)
<u>23 ml 1X PBS</u>	
25 ml total	

4. Rinse the cells twice with 1X PBS.

5. Add the Staining Solution to the cells and incubate at 37°C for 30 minutes to 2 hours.
6. Check the cells under a microscope. Calculate the percent of cells transfected with  $\beta$ -galactosidase.

$$\frac{\text{Total No. of blue cells}}{\text{Total No. of cells}} \times 100 = \% \text{ transfection}$$

## References

---

1. Capecchi, M.R. (1980) *Cell* 22: 479-488.
2. Graham, F.L. and van der Eb, A.J. (1973) *Virology* 52: 456-467.
3. Felgner, P.L. *et al.* (1987) *Proc. Natl. Acad. Sci. USA* 84: 7413-7417.
4. Cepko, C.L. *et al.* (1984) *Cell* 37: 1053-1062.
5. Neumann, E. *et al.* (1982) *EMBO J.* 7: 841-845.
6. Schwarze, S.L. *et al.* (2000) *Trends Cell Biol.* 10: 290-295.
7. Morris, M. (1999) *et al. J. Biol. Chem.* 274 (35): 24941-24946.
8. Morris, M. *et al.* (2001) *Nature Biotech.* 19: 1173-1176.

## Troubleshooting Guide

PROBLEM	POSSIBLE CAUSE	RECOMMENDATION
Low Transfection Efficiency	Cell density	Use adherent cells at a confluency of 40-50%.
	Chariot or macromolecule dilutions were performed in larger volumes than recommended	Dilute the macromolecule into the final volume of PBS as recommended. Dilute Chariot in the volume of sterile water recommended. It is necessary to make the Chariot macromolecule complex in a concentrated solution. For multiple transfections, do not use a master mix that exceeds the volume required for 4 transfections (6-well plate size) per tube, as this may cause aggregation.
	Chariot and macromolecule were not complexed for 30 minutes	Combine the Chariot dilution and the macromolecule dilution together and allow the complex to form at room temperature for 30 minutes.
	Improper storage of Chariot	Chariot may form aggregates over time if not stored properly. You may need to sonicate the Chariot dilution just before complex formation. We recommend placing the Chariot dilution in a water bath and sonicating the water at close proximity with an ultrasonic processor. At Active Motif, we routinely use Sonics Vibra-cell™ for 1 minute at an amplitude of 30 and a 1-second pulse.
	Medium that contained serum was used to dilute the Chariot-macromolecule complex	Repeat the transfection using serum-free medium to dilute the Chariot-macromolecule complex to the Final Transfection Volume.
	Exposure of Chariot-macromolecule complex to cells was insufficient	Leave the Chariot-macromolecule complex on the cells for a minimum of two hours.
	Using a cell line other than those tested with Chariot	Try the transfection in a cell line proven to work with Chariot. Please visit our website for complete information on cell lines that have been tested
	A suboptimal amount of macromolecule was used for the transfection	You may need to titrate the amount of macromolecule (protein, peptide or antibody) used in the transfection.
A suboptimal amount of Chariot was used for the transfection	You may need to titrate the amount of Chariot used in the transfection. Chariot interacts via hydrophobic interactions. Each peptide or protein will have a different hydrophobicity and may require a different amount of Chariot. Depending on the sensitivity required for antibody detection it may be necessary to alter the amount of Chariot used. Detection of fluorescently labeled antibodies in cells requires a large amount of antibody to be delivered. This may mean increasing the amount of Chariot used.	
Diluting Chariot into DMSO may improve antibody delivery	Diluting Chariot into DMSO has improved the delivery of some antibodies. If delivery of your antibody is low, try diluting Chariot into 60-80% DMSO instead of water. After the Chariot dilution is combined with the antibody dilution, the complex mixture will have a final concentration of 30-40% DMSO. This may need to be lowered for certain cell lines. However, a final concentration of complex mixture below 10-20% DMSO has not been shown to improve antibody delivery.	

PROBLEM	POSSIBLE CAUSE	RECOMMENDATION
Protein delivery is localized to a few regions of the cell	Diluting Chariot into DMSO may make protein delivery more uniform	With some proteins that we have tested, delivery of the protein seems to be localized to specific regions in the cell. Diluting Chariot into DMSO instead of water has made the delivery of some of these proteins more uniform. Dilute the Chariot into 40-60% DMSO. After the Chariot dilution is combined with the protein dilution, the complex mixture will have a final concentration of 20-30% DMSO.
Aggregate Formation	Improper storage of Chariot	Chariot may form aggregates over time if not stored properly. You may need to sonicate the Chariot dilution just before complex formation. We recommend placing the Chariot dilution in a water bath and sonicating the waterbath with an ultrasonic processor, such as a Sonics Vibra-cell™, for 1 minute at an amplitude of 30 and a 1-second pulse.
	Excess Chariot used	Decrease the amount of Chariot used
Signs of cytotoxicity. <b>Note:</b> Chariot is non-cytotoxic to cells at the recommended concentrations.	Transfected macromolecule may be cytotoxic	Transfect the macromolecule at a lower concentration. Compare cells only, cells with Chariot alone and cells with macromolecule alone. Chariot should not affect cell viability. Transfect the β-galactosidase control protein.
	Culture may be contaminated with mycoplasma	Treat the cells to eliminate mycoplasma.
	Cells may not be healthy	Check the incubator (CO <sub>2</sub> and temperature levels). Check the media.
	Too much DMSO present	Reduce the final concentration of DMSO to a final concentration below 40%.

## Related Products

Fluorescent Dyes	Units	Catalog No.
Chromo™ 494 Carboxylic Acid	1 mg	15110
	5 mg	16110
Chromo™ 494 NHS Ester	1 mg	15111
	5 mg	16111
Chromo™ 494 Biotin	1 mg	15112
	5 mg	16112
Chromo™ 494 Streptavidin	1 mg	15113
	5 mg	16113
Chromo™ 546 Carboxylic Acid	1 mg	15210
	5 mg	16210
Chromo™ 546 NHS Ester	1 mg	15211
	5 mg	16211
Chromo™ 546 Biotin	1 mg	15212
	5 mg	16212
Chromo™ 546 Streptavidin	1 mg	15213
	5 mg	16213
Chromo™ 642 Carboxylic Acid	1 mg	15310
	5 mg	16310
Chromo™ 642 NHS Ester	1 mg	15311
	5 mg	16311
Chromo™ 642 Biotin	1 mg	15312
	5 mg	16312
Chromo™ 642 Streptavidin	1 mg	15313
	5 mg	16313

## 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](mailto:tech_service@activemotif.com)

### Active Motif Europe

104 Avenue Franklin Roosevelt

B-1330 Rixensart, 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](mailto: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](mailto:japantech@activemotif.com)

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, *Motivations* 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!**