

LigandLink™

Universal Labeling Technology

(version A1)

Product	Catalog No.
LigandLink™ pLL-1 Kit	34001
Premade LigandLink™ Kits	34002-34006
LigandLink™ Fluorescein Label	34101
LigandLink™ Hexachlorofluorescein Label	34104

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

Background	
Overview . . .	1
Introduction	2
LigandLink pLL-1 Vector Map	3
Sequence of pLL-1 MCS	4
LigandLink pLL-1 NLS Vector Map	5
Labeling of nuclear localized eDHFR using LigandLink	6
Kit Components and Storage	7
Additional Materials Required	8
Component Preparation and Recommendations	9
LigandLink Cell Labeling Protocols	
Mammalian Cell Transfection	10
Cell Staining with LigandLink Label	11
References	11
Appendix	
Section A. Troubleshooting	12
Section B. Related Products	13
Technical Services	14

Overview

Active Motif's LigandLink™ Universal Labeling Kit enables specific and flexible labeling of proteins in living cells. In the LigandLink method, the gene of interest is cloned in frame with the gene for *E. coli* dihydrofolate reductase (eDHFR) in the pLL-1 vector. The vector is then transfected into mammalian cells and used to express fusion protein. After transfection, the protein of interest can be labeled simply by adding the LigandLink Label of choice to the cell medium (Figure 1). Depending on the cell type and the characteristics of the label, cells can be imaged in as little as ten minutes.

product	format	catalog no.
LigandLink™ pLL-1 Kit	1 kit	34001
Premade LigandLink™ Kit	1 kit	34002-34006
LigandLink™ Fluorescein Label	300 rxns	34101
LigandLink™ Hexachlorofluorescein Label	300 rxns	34104

Each LigandLink Kit contains 20 µg of vector, 20 µg of LigandLink pLL-1 NLS positive control, 100 reactions of LigandLink Fluorescein Label and one vial of DMF + acetic acid. Each reaction of LigandLink Label is sufficient to label a single well of a 96-well plate.

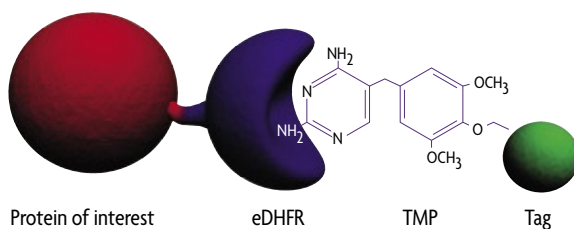


Figure 1: Specific protein labeling using LigandLink.

The gene of interest is cloned into pLL-1 in frame with the vector's *E. coli* dihydrofolate reductase (eDHFR) gene. After transfection into cells, the protein of interest is expressed as a fusion to eDHFR. Addition of cell-permeable LigandLink Label to the medium results in rapid, specific binding of the label by the fusion protein.

Introduction

Use of green fluorescent protein (GFP) has become commonplace in modern biology. However, biologically fluorescent proteins (FP) are limited by their inherent properties. For example, FPs have a relatively low quantum yield, and it is difficult to engineer their spectral properties to suit specific applications. Fluorescent dyes are a potential alternative to biological FPs as they are available in a broad variety of formats and can be engineered easily to ensure desired spectral properties. However, because fluorescent dyes are synthetic molecules it has not been possible to use them as a general tool for labeling specific proteins within a cell.

Active Motif's LigandLink Universal Labeling technology overcomes these limitations by providing a small ligand that can carry a variety of functional tags, including fluorescent dyes. Expression of your protein of interest as a fusion enables you to specifically label the protein of interest *in vivo* simply by adding one of the LigandLink labels to the medium*. To change the properties of your tag all you need to do is add a different LigandLink label. Thus LigandLink enables you to create a single protein fusion that can be labeled with a variety of tags, depending on the needs of your experiment.

The eDHFR protein was chosen as a fusion protein because it is a relatively small, monomeric protein (18 kDa vs 27 kDa for GFP) that has been shown to have a high affinity for the ligand trimethoprim (TMP). TMP binds with a high specificity to the *E. coli* form of DHFR ($K_1 = -1$ nm), and a substantially lower affinity for endogenous DHFR ($K_1 = -4$ μ m). This is because TMP is an antibiotic that was designed to specifically inhibit the bacterial enzymes responsible for the production of folic acid while not interacting with mammalian proteins. As there is minimal binding of LigandLink Labels to non-tagged mammalian proteins, this results in extremely low background. Moreover, TMP can be derivatized to carry a number of tags without substantially altering its affinity and specificity for eDHFR.

The LigandLink Universal Labeling technology uses eDHFR and a number of TMP derivatives as a ligand-receptor pair to provide a variety of functionalities¹. Cell-permeable LigandLink Labels are available with Fluorescein and Hexachlorofluorescein fluorescent dyes.

* Patent pending.

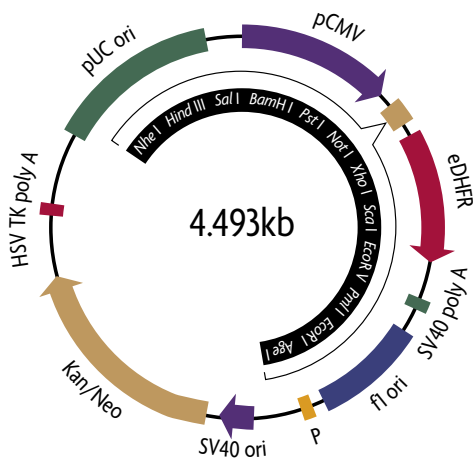


Figure 2: The LigandLink pLL-1 vector.

The LigandLink vector, pLL-1, was designed for ease of use. It features a CMV promoter for high level expression of eDHFR fusion proteins, with Neomycin for selection of stable cell lines. The multiple cloning site (MCS) was designed to facilitate cloning, whatever method you use. In addition to many popular restriction sites, the MCS includes three blunt-cutting restriction enzymes towards the 3' end, each in a different reading frame with the eDHFR gene. This enables a number of PCR and restriction enzyme cloning strategies.

pLL-1 Vector Sequence Reference Points

Human cytomegalovirus (CMV) immediate early promoter	1-589
MCS	591-662
eDHFR	670-1170
SV40 early mRNA polyadenylation signal	1312-1345
f1 single-strand DNA origin	1409-1864
Bacterial promoter for expression of Kan ^r gene	1926-1954
SV40 origin of replication	2205-2340
Kanamycin/neomycin resistance gene	2389-3183
Herpes simplex virus (HSV) thymidine kinase (TK) polyadenylation signal	3420-3437
pUC plasmid replication origin	3768-4411
pLL-1 FWD: 5' -ATGTCGTAACAACCTCCGCC-3'	493-512
pLL-1 REV: 5' -CAGGTTCCACGGCATGGCGTT-3'	742-762

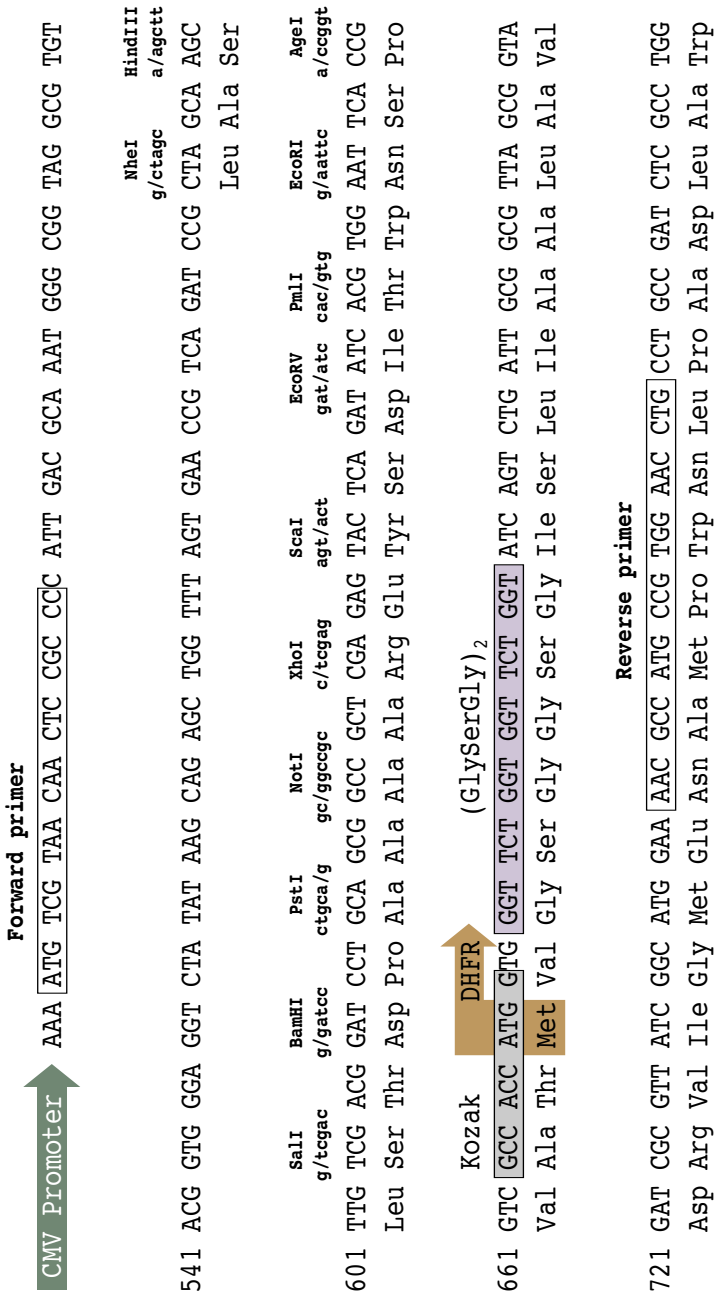


Figure 3: The sequence of the pLL-1 vector illustrating the cloning and primer hybridization sites.

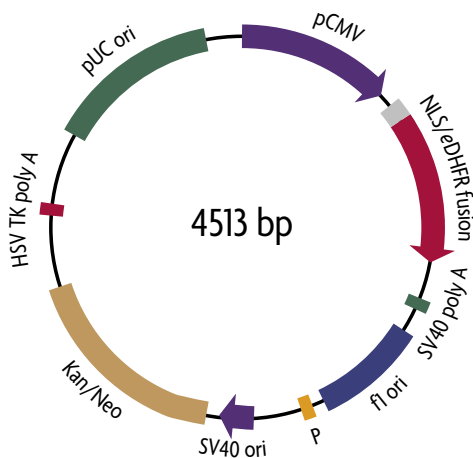


Figure 4: The LigandLink pLL-1 NLS vector.

To generate pLL-1 NLS, three copies of the canonical simian virus 40 T-antigen nuclear localization sequence (DPKKRKRKV) were fused to the N-terminus of eDHFR in pLL-1 by preparing a double stranded oligonucleotide and ligating the NLS sequence into the *Nhe* I and *Age* I sites. This vector can be used as a positive control for nuclear staining as the NLS/eDHFR protein will translocate to the nucleus.

pLL-1 NLS Vector Sequence Reference Points

Human cytomegalovirus (CMV) immediate early promoter	1-589
Nuclear localization site (NLS) coding sequence:	603-677
eDHFR	690-1190
SV40 early mRNA polyadenylation signal	1332-1365
f1 single-strand DNA origin	1429-1864
Bacterial promoter for expression of Kan ^r gene	1946-1974
SV40 origin of replication	2205-2340
Kanamycin/neomycin resistance gene	2409-3203
Herpes simplex virus (HSV) thymidine kinase (TK) polyadenylation signal	3440-3457
pUC plasmid replication origin	3788-4431
pLL-1 FWD: 5´-ATGTCGTAACAACCTCCGCCC-3´	493-512
pLL-1 REV: 5´-CAGGTTCCACGGCATGGCGTT-3´	762-782

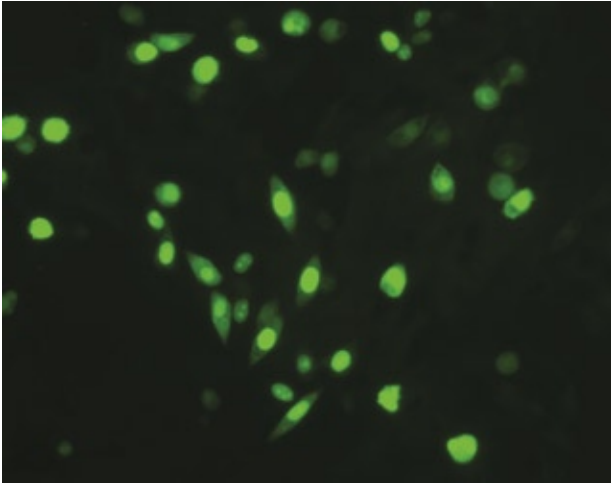


Figure 5: Labeling of nuclear localized eDHFR by LigandLink Fluorescein.

A nuclear localization sequence was cloned into pLL-1 and transfected into CHO cells. Twenty-four hours post transfection, 5 μ M LigandLink Fluorescein was added to the cells for 2 hours at 37°C. Following wash steps to remove unbound label, the above image was taken.

Kit Components and Storage

Kit components can be stored at -20°C prior to first use. Then, we recommend storing each component at the temperature indicated in the table below.

LigandLink pLL-1 Kit	Quantity	Storage / Stability
pLL-1 Vector	20 µg	-20°C for 6 months
pLL-1 NLS (positive control vector for nuclear translocation)	20 µg	-20°C for 6 months
LigandLink Fluorescein Label	100 rxns	-20°C for 6 months
DMF + acetic acid	50 µl	4°C for 6 months

Premade LigandLink Kits	Quantity	Storage / Stability
pLL-1 eDHFR Fusion	20 µg	-20°C for 6 months
pLL-1 NLS (positive control vector for nuclear translocation)	20 µg	-20°C for 6 months
LigandLink Fluorescein Label	100 rxns	-20°C for 6 months
DMF + acetic acid	50 µl	4°C for 6 months

LigandLink Labels	Quantity	Storage / Stability
LigandLink Fluorescein Label*	300 rxns	-20°C for 6 months
LigandLink Hexachlorofluorescein Label	300 rxns	-20°C for 6 months

* Each order of LigandLink Label is supplied with DMF + acetic acid (150 µl). This may be stored at 4°C for 6 months. Each reaction of LigandLink Label is sufficient to label a single well of a 96-well plate.

The LigandLink Universal Labeling Kit is for research use only. Not for use in diagnostic procedures.

Additional Materials Required

Multi-channel pipettor
Multi-channel pipettor reservoirs
Chambered covered glass slide or 96-well plate (or alternate seeding plate)
Transfection Reagent
Endotoxin-free (transfection-grade) plasmid DNA
Fetal bovine serum (FBS)
Serum-free cell culture medium
Phosphate buffered saline (PBS, pH 7.4) containing Ca^{2+} and Mg^{2+}
Glucose
Sterile distilled water (dH_2O)
Fluorescent microscope equipped with standard FITC filters
37°C cell culture incubator

Protocols

Component Preparation and Recommendations

Prior to starting the assay please prepare the following:

Preparation of Plasmid DNA

Stock Solution: The pLL-1, pLL-1 NLS control, and all premade pLL-1 vectors are supplied lyophilized as 20 µg aliquots. Prepare the DNA Stock Solution by resuspending the lyophilized DNA in 20 µl of sterile dH₂O in the provided vial. Stock concentration is 1 µg/µl. This Stock Solution can be stored at -20°C for 6 months.

Working Solution: Prepare the amount of plasmid DNA required for the assay by diluting the plasmid DNA in sterile dH₂O. For example for an 8 well chamber slide mix 1.5 µl of stock DNA with 23.5 µl of sterile dH₂O to generate a 0.06 µg/µl Working Solution.

Resuspension of LigandLink Fluorescent Label

The LigandLink Fluorescent Label is supplied lyophilized. The quantity of label supplied is 30 nmoles*. Prepare the LigandLink Fluorescent Label Stock Solution by resuspending the lyophilized label in 30 µl of DMF + acetic acid in the provided amber vial. This generates a 1 mM stock solution. This Stock Solution can be stored at -20°C for 6 months. **Note:** Fluorescent dyes are light sensitive. Avoid exposing the cells to light during the ligand labeling and wash steps.

* For the 300-reaction aliquots of LigandLink Labels, three vials each with 30 nmoles of lyophilized label are provided.

Mammalian Cell Transfection

These conditions are recommended as guidelines only. The protocol below has been optimized for transfection of CHO-K1 cells (ATCC Cat. No. CCL-61) in an Lab Tek® 8-well chambered covered glass slide (Nalge Nunc Cat. No. 177402) using FuGENE® 6 Transfection Reagent (Roche Cat. No. 11 814 443 001). Please note that other transfection reagents may be used. You should empirically optimize the cell culture protocol, transfection conditions, ligand concentration and labeling protocol for your experimental system.

1. In an 8-well chambered covered glass slide, seed 1.6×10^4 cells per well in 250 μ l of complete growth medium.
2. Incubate the cells at 37°C in a humidified atmosphere containing 5% CO₂ until the cells are 90-95% confluent. For CHO-K1 cells this requires 16-20 hours.
3. In a microcentrifuge tube mix 88 μ l of a serum-free culture media without additives or supplements with 4.4 μ l FuGENE® 6 Transfection Reagent. **Note:** Do not touch the plastic walls of the tubes.
4. Tap the tube gently. Incubate at room temperature for 5 minutes.
5. Add 22 μ l of the working DNA solution (see Preparation of Reagents above).
6. Tap the tube gently. Incubate at room temperature for 15 minutes.
7. Tap the tube gently again. Overlay 10 μ l onto each well.
8. Incubate cells at 37°C in a humidified atmosphere containing 5% CO₂ overnight.

Cell Staining with LigandLink Label

This staining protocol is suitable for both LigandLink Fluorescein Label and LigandLink Hexachlorofluorescein Label.

Note 1: After removing the LigandLink Label from -20°C, we recommend placing it at room temperature for 5 minutes before use. This minimizes moisture from entering the vial upon opening.

Note 2: Fluorescent dyes are light sensitive. Avoid exposing the cells to light during the labeling and wash steps.

Note 3: Only prepare the amount of diluted label required for the assay and use only freshly prepared diluted LigandLink Label.

Note 4: The concentration of LigandLink Label, the time of labeling, the wash conditions and the microscope exposure need to be optimized for each cell line.

Note 5: LigandLink Labels can be added directly to serum-containing medium.

Note 6: Use a standard FITC filter set to detect proteins labeled with LigandLink Fluorescein Label. Excitation of 488 nm and emission between 500 and 550 nm. The optimal excitation and emission wavelengths for LigandLink Hexachlorofluorescein Label are 535 and 565 nm, respectively.

Note 7: Due to the acidity of the resuspended label, a slight color change may be observed in media that contains phenol red. This does not affect the results.

1. Dilute the LigandLink Label stock solution 1-5 μM in cell medium.
2. Add 200 μl of diluted label to each well of the 8-well chambered covered glass slide.
3. Incubate 2 hours at 37°C in a humidified atmosphere containing 5% CO_2 .
4. Wash cells twice with 200 μl of PBS.
5. Add 200 μl of PBS supplemented with 10 mM glucose to each well of the 8-well chambered covered glass slide. We recommend that you use a PBS that contains both Ca^{2+} (-0.9 mM) and Mg^{2+} (-0.49 mM) to maintain healthy cells.
6. Transfer the chambered cover glass slide to a microscope and capture images.

References

1. Miller L.W. *et al.* (2005) *Nature Methods* 2(4): 255-257.

Appendix

Section A. Troubleshooting Guide

PROBLEM	POSSIBLE CAUSE	RECOMMENDATION
No signal or weak signal	Fusion protein not expressed or expressed only at low levels	Check the reading frame of your construct by sequence analysis. Optimize transfection conditions and use high-quality, endotoxin-free DNA. Culture cells for a longer period of time before labeling to ensure that you have adequate protein expression. Protein expression can be improved by optimizing the health of the cells. Increase the seeding density or time of culture to allow cells to proliferate and adhere more tightly.
	Fusion protein stop codon not removed prior to cloning into pLL-1	Because the pLL-1 vector contains an MCS that is upstream of the eDHFR coding sequence, make sure the stop codon is removed from the protein of interest before cloning to ensure that the gene of interest will be in frame with the downstream eDHFR coding sequence (which contains its own stop codon).
	Insufficient LigandLink Label	Optimize cell-labeling protocols. Increase cell labeling time. Optimize the ligand concentration. Increase the exposure time while viewing cells under the microscope.
	Loss of labeling activity	Store the LigandLink Labels at -20°C and protect them from light so that they do not lose labeling activity. Use freshly prepared labeling solution to label your cells.
	Incorrect filter set	Ensure that you are using the correct filter set for viewing. Use a standard FITC filter set to detect proteins labeled with the LigandLink Fluorescein Label. Excitation of 488 nm and emission between 500 and 550 nm. The optimal excitation and emission wavelengths for LigandLink Hexachlorofluorescein Label are 535 and 565 nm respectively.
	Ligand photobleaching	To prevent LigandLink Labels from photobleaching, analyze the fluorescent signal for only a short period of time.
High background fluorescence	Inadequate washing	Increase time for washing unbound ligand. Reduce label concentration. Reduce labeling time. Reduce exposure time while viewing cells under the microscope.
Cells detach from surface		When imaging live cells, handle cells carefully to ensure that they stay attached to the surface.
Altered cell morphology		Label cells in the dark so they are not exposed to intense light during labeling.

Section B. Related Products

TransAM™ Kits	Unit	Catalog No.	
TransAM™ NFκB Family	2 x 96 rxns	43296	
TransAM™ Flexi NFκB Family	2 x 96 rxns	43298	
TransAM™ NFκB p50	1 x 96 rxns	41096	
	5 x 96 rxns	41596	
TransAM™ NFκB p50 Chemi	1 x 96 rxns	41097	
	5 x 96 rxns	41597	
TransAM™ Flexi NFκB p50	1 x 96 rxns	41098	
TransAM™ NFκB p65	1 x 96 rxns	40096	
	5 x 96 rxns	40596	
TransAM™ NFκB p65 Chemi	1 x 96 rxns	40097	
	5 x 96 rxns	40597	
TransAM™ Flexi NFκB p65	1 x 96 rxns	40098	
TransAM™ STAT Family	2 x 96 rxns	42296	
TransAM™ STAT3	1 x 96 rxns	45196	
	5 x 96 rxns	45696	
Cell-based ELISAs	Unit	Colorimetric Kit Catalog No.	Chemi Kit Catalog No.
FACE™ AKT	1 x 96 rxns	48120	48220
	5 x 96 rxns	48620	48720
FACE™ NFκB p65 Profiler	3 x 96 rxns	48300	48400
FACE™ STAT2	1 x 96 rxns	48310	48410
	5 x 96 rxns	48810	48910
FACE™ STAT4	1 x 96 rxns	48320	48420
	5 x 96 rxns	48820	48920
FACE™ STAT6	1 x 96 rxns	48330	48430
	5 x 96 rxns	48830	48930
Competent Cells	Unit	Catalog No.	
RapidTrans™ TAMI-F´ Competent <i>E. coli</i>	1 x 96 rxns	10096	
	5 x 96 rxns	10596	
RapidTrans™ TAMI Competent <i>E. coli</i>	1 x 96 rxns	11096	
	5 x 96 rxns	11596	
RapidTrans™ TAMI λ pir+ Competent <i>E. coli</i>	1 x 96 rxns	11097	
	5 x 96 rxns	11597	
RapidTrans™ TAP Competent <i>E. coli</i>	1 x 96 rxns	11098	
	5 x 96 rxns	11598	
RapidTrans™ TAP-F´ Competent <i>E. coli</i>	1 x 96 rxns	10098	
	5 x 96 rxns	10598	
Protein Transfection	Unit	Catalog No.	
Chariot™	25 rxns	30025	
	100 rxns	30100	
Fluorescent Dyes	Unit	Catalog No.	
Chromo™ 494 Carboxylic Acid	1 mg	15110	
	5 mg	16110	
Chromo™ 494 NHS-Ester	1 mg	15111	
	5 mg	16111	
Chromo™ 546 Carboxylic Acid	1 mg	15210	
	5 mg	16210	
Chromo™ 546 NHS-Ester	1 mg	15211	
	5 mg	16211	
Chromo™ 642 Carboxylic Acid	1 mg	15310	
	5 mg	16310	
Chromo™ 642 NHS-Ester	1 mg	15311	
	5 mg	16311	

Technical Services

If you need assistance at any time, please call the Active Motif Technical Service Department 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

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

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

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!