

Click Chemistry Reagents

Table 1. Contents and storage information.

Material	Amount	Storage	Stability
Alkyne or azide-containing click chemistry reagent	Varies, see product label	≤-20°CDesiccateProtect from light	When stored as directed the product is stable for 6–12 months.
Approximate fluorescence excitation/emiss	ion maxima: See Table 2.		

Introduction

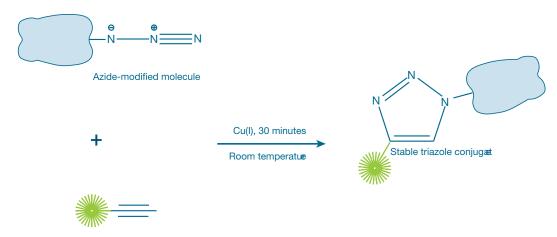
Click chemistry describes a class of chemical reactions that use bio-orthogonal or biologically unique moities to label and detect a molecule of interest using a two-step procedure.¹⁻⁴ The two-step click reaction involves a copper-catalyzed triazole formation from an azide and an alkyne (Figure 1). The azide and alkyne moieties can be used interchangeably; either one can be used to tag the molecule of interest, while the other is used for subsequent detection. The azides and alkynes are biologically unique, inert, stable, and extremely small (Figure 2).

Click chemistry can be used when methods such as direct labeling or the use of antibodies are not applicable or efficient. The click chemistry label is small enough that tagged molecules (e.g., nucleotides⁵, sugars⁶, and amino acids⁷) are acceptable substrates for the enzymes that assemble these building blocks into biopolymers. The small size of click detection molecules allows them to easily penetrate complex samples, including intact, supercoiled DNA, with only mild permeabilization required.

The characteristics of click reactions include:

- Efficiency—the reaction between the detection moieties is complete in less than 1 hour and does not require extreme temperatures or solvents.
- **Stability**—the reaction product contains an irreversible, covalent bond.
- Biologically inert—the components of the reaction do not undergo any side reactions.
- **Specificity**—the reaction between the label and detection tag is selective and specific.
- Applicability to biological samples—the click chemistry-labeled molecules can be applied to complex biological samples and easily detected with high sensitivity and low background, unlike traditional chemical reactions that use succinimidyl esters or maleimides that target amines and sulfhydryls, which are not unique functional groups.

A wide variety of azide- or alkyne-containing dyes, haptens, and biomolecules for use in click reactions are available from Invitrogen (Tables 2-4). A general protocol for the coppercatalyzed click reaction between an azide and an alkyne is described below.



Fluorophore-, or hapten-alkyne

Figure 1. Click azide/alkyne reaction. The azide and alkyne moieties are interchangeable, whereupon the molecule can be labeled with an alkyne and reacted with a fluorophore or hapten-azide.

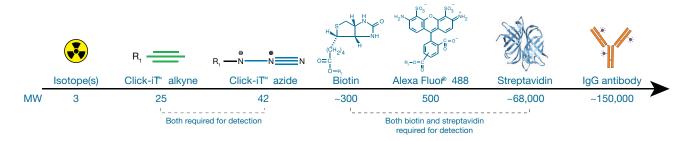


Figure 2. Relative size of detection molecules commonly used in cellular analysis.

Table 2. Azide- or alkyne-modified fluorophores and haptens.

Label	Ex/Em*	Azide or alkyne	Cat. no.	Use
Alexa Fluor® 488	495/519	azide	A10266	Fluorescent due or hanten
Alexa Fluor® 488	495/519	alkyne	A10267	Fluorescent dye or hapten
Alexa Fluor® 555	FF /F /F	azide	A20012	
Alexa Fluor® 555	555/565	alkyne	A20013	
Alexa Fluor® 594	500/617	azide	A10270	Fluores and due
Alexa Fluor® 594	590/617	alkyne	A10275	Fluorescent dye
Alexa Fluor® 647	650/655	azide	A10277	
Alexa Fluor® 647	650/655	alkyne	A10278	
Distin	Not	azide	B10184	11
Biotin	applicable	alkyne	B10185	Hapten
Out 11.00 City 11.00	406/524	azide	O10180	
Oregon Green® 488	496/524	alkyne	O10181	
Takan and all design and and and	FFF /F00	azide	T10182	Fluorescent dye or hapten
Tetramethylrhodamine	555/580	alkyne	T10183	
*Fluorescence excitation	and emission m	axima in nm.		

Table 3. Azide- or alkyne-modified biomolecules.

Compound	Cat. no.	Azide or alkyne	Application
EdU (5'-ethynyl-2'-deoxyuridine)	A10044, E10187	alkyne	For nascent DNA synthesis
Click-iT® AHA (L-azidohomoalanine)	C10102	azide	Nascent protein synthesis
Click-iT® HPG (L-homopropargylglycine)	C10186	alkyne	Nascent protein synthesis
Click-iT® farnesyl alcohol, azide	C10248	azide	Farnesylated proteins
Click-iT® geranylgeranyl alcohol, azide	C10249	azide	Geranylgeranylated proteins
Click-iT® fucose alkyne (tetraacetylfucose alkyne)	C10264	alkyne	Fucosylated glycans
Click-iT® palmitic acid, azide	C10265	azide	Palmitoylated proteins
Click-iT® myristic acid, azide	C10268	azide	Myristoylated proteins
Click-iT® GalNAz (tetraacetylated N-azidoacetylgalactosamine)	C33365	azide	O-Linked glycoproteins
Click-iT® ManNAz (tetraacetylated N-azidoacetyl-D-mannosamine)	C33366	azide	Sialic acid-modified glycoproteins
Click-iT® GlcNAz (tetraacetylated N-azidoacetylglucosamine)	C33367	azide	O-GlcNAz-modified glycoproteins

 Table 4. Reactive azides and alkynes.

Reactivity	Reactive moiety	Azide or alkyne	Cat. no.
Duineauraninaa	Consissional deal and an	azide	A10280
Primary amines	Succinimidyl ester	alkyne	A10279
Thirds	la da casta mila	azide	I10188
Thiols	lodoacetamide	alkyne	I10189

Before Starting

Azide- and Alkyne-modified-Molecules

The key components to any click reaction are an azide labeled molecule and an alkyne labeled molecule.

Solvent

- Most alkyne- and azide-modified fluorophores, haptens, and reactive probes are hydrophobic molecules. We recommend that you dissolve these molecules in high-quality, anhydrous dimethylformamide (DMF) or dimethylsulfoxide (DMSO).
- For azide- and alkyne-modified biomolecules (Table 3), refer to the specific protocol supplied with the products.

Click Reaction Conditions

Use of Copper as a Catalyst

In addition to the azide and alkyne labeled molecules, copper (I) is required to catalyze the reaction. We recommend using copper (II) sulfate in the presence of a reductant such as ascorbic acid to generate copper (I). The use of copper (I) directly is less favored due to the ease with which it is oxidized to the non-catalytic copper (II) species. The preferred method is the reduction of copper (II) sulfate *in situ* to obtain copper (I).

pH and Temperature

The click reaction is highly efficient and extremely tolerant of a wide variety of conditions. The click reaction occurs at pH values ranging from 3 to 12, at room temperature, generally in less than one hour. Reaction rates have been found to increase slightly at lower pH levels where copper (I) is more soluble, and therefore more readily available for catalysis. Although the reaction is not light sensitive, protect the reaction from light for light-sensitive fluorophores.

Click-iT® Reaction Buffer Kits

For convenience, Invitrogen offers Click-iT* Reaction Buffer Kits for protein or cell samples labeled with an azide- or alkyne-tagged biomolecule. The Click-iT° Cell Reaction Buffer Kit (Cat. no. C10269) includes sufficient reagents to perform 50 reactions based on a 0.5 mL reaction volume for subsequent analyses by flow cytometry, fluorescence microscopy, or high content screening (HCS). The Click-iT® Protein Reaction Buffer Kit (Cat. no. C10276) includes everything required to perform the click reaction of proteins for subsequent standard protein biochemical analyses (e.g., western blots, mass spectrometry).

References

1. ChemBioChem 4, 1147 (2003); 2. J Am Chem Soc 125, 3192 (2003); 3. Angew Chem Int Ed Engl 41, 2596 (2002); 4. Angew Chem Int Ed Engl 40, 2004 (2001); 5. Proc Natl Acad Sci 105, 2415 (2008); 6. J Am Chem Soc 130, 11576 (2008); 7. Proc Natl Acad Sci 103, 9482 (2006).

Product List Current prices may be obtained from our website or from our Customer Service Department.

Cat. no.Product NameUnit SizeA10266Alexa Fluor* 488 azide (Alexa Fluor* 488 5-carboxamido-(6-azidohexanyl), bis(triethylammonium salt)).0.5 mgA10267Alexa Fluor* 488 alkyne (Alexa Fluor* 488 5-carboxamido-(propargyl), bis(triethylammonium salt)).0.5 mgA20012Alexa Fluor* 555 azide, triethylammonium salt.0.5 mgA20013Alexa Fluor* 555 alkyne, triethylammonium salt.0.5 mgA10270Alexa Fluor* 594 azide (Alexa Fluor* 594 carboxamido-(6-azidohexanyl), bis(triethylammonium salt)).0.5 mgA10275Alexa Fluor* 594 alkyne (Alexa Fluor* 594 carboxamido-(5-(and 6-)propargyl), bis(triethylammonium salt)).0.5 mgA10277Alexa Fluor* 647 alkyne, triethylammonium salt.0.5 mgA10278Alexa Fluor* 647 alkyne, triethylammonium salt.0.5 mgA10279alkyne, succinimidyl ester (3-propargyloxypropanoic acid, succinimidyl ester).1 mgB10184biotin azide.1 mgB10185biotin alkyne.1 mgB10188biodacetamide azide.1 mgB10189iodoacetamide alkyne.1 mgB10180Oregon Green* 488 alkyne *6-isomer*0.5 mgD10181Oregon Green* 488 alkyne *6-isomer*0.5 mgT10182tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer*0.5 mgT10183tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer*0.5 mgA10044EdU (5-ethynyl-2'-deoxyuridine)50 mgC10102Click-iT* AHA (L-azidohomoalanine) *for nascent protein synthesis* <t< th=""></t<>
A10267 Alexa Fluor® 488 alkyne (Alexa Fluor® 488 5-carboxamido-(propargyl), bis(triethylammonium salt)). 0.5 mg A20012 Alexa Fluor® 555 azide, triethylammonium salt . 0.5 mg A20013 Alexa Fluor® 555 alkyne, triethylammonium salt . 0.5 mg A10270 Alexa Fluor® 594 azide (Alexa Fluor® 594 carboxamido-(6-azidohexanyl), bis(triethylammonium salt)). 0.5 mg A10275 Alexa Fluor® 647 azide, triethylammonium salt . 0.5 mg A10276 Alexa Fluor® 647 azide, triethylammonium salt . 0.5 mg A10277 Alexa Fluor® 647 azide, triethylammonium salt . 0.5 mg A10278 Alexa Fluor® 647 azide, triethylammonium salt . 0.5 mg A10279 alkyne, succinimidyl ester (3-propargyloxypropanoic acid, succinimidyl ester). 1 mg B10184 biotin azide . 1 mg B10185 biotin alkyne. 1 mg B10188 iodoacetamide azide . 1 mg B10189 iodoacetamide azide . 1 mg B10180 Oregon Green® 488 azide (Oregon Green® 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt) . 0.5 mg C10181 Oregon Green® 488 alkyne %6-isomer® . 0.5 mg Related Products A10044 EdU (5-ethynyl-2'-deoxyuridine) . 50 mg C10102 Click-iT® AHA (L-azidohomoalanine) *for nascent protein synthesis® . 5 mg
A20012Alexa Fluor® 555 azide, triethylammonium salt0.5 mgA20013Alexa Fluor® 555 alkyne, triethylammonium salt0.5 mgA10270Alexa Fluor® 594 azide (Alexa Fluor® 594 carboxamido-(6-azidohexanyl), bis(triethylammonium salt))0.5 mgA10275Alexa Fluor® 594 alkyne (Alexa Fluor® 594 carboxamido-(5-(and 6-)propargyl), bis(triethylammonium salt))0.5 mgA10277Alexa Fluor® 647 azide, triethylammonium salt0.5 mgA10278Alexa Fluor® 647 alkyne, triethylammonium salt0.5 mgA10279alkyne, succinimidyl ester (3-propargyloxypropanoic acid, succinimidyl ester)1 mgA10280azido (PEO)₄ propionic acid, succinimidyl ester (3-(azidotetra(ethyleneoxy))propionic acid, succinimidyl ester)1 mgB10184biotin azide1 mgB10185biotin alkyne1 mgI10188iodoacetamide alkyne1 mgI10189iodoacetamide alkyne1 mg010180Oregon Green® 488 azide (Oregon Green® 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt)0.5 mg010181Oregon Green® 488 alkyne ®6-isomer®0.5 mgT10182tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) ®5-isomer®0.5 mgRelated Products50 mgA10044EdU (5-ethynyl-2'-deoxyuridine)50 mgClick-iT® AHA (L-azidohomoalanine) *for nascent protein synthesis®5 mg
A20013 Alexa Fluor* 555 alkyne, triethylammonium salt. 0.5 mg A10270 Alexa Fluor* 594 azide (Alexa Fluor* 594 carboxamido-(6-azidohexanyl), bis(triethylammonium salt)) 0.5 mg A10275 Alexa Fluor* 594 alkyne (Alexa Fluor* 594 carboxamido-(5-(and 6-)propargyl), bis(triethylammonium salt)) 0.5 mg A10277 Alexa Fluor* 647 azide, triethylammonium salt 0.5 mg A10278 Alexa Fluor* 647 alkyne, triethylammonium salt 0.5 mg A10279 alkyne, succinimidyl ester (3-propargyloxypropanoic acid, succinimidyl ester) 1 mg A10280 azido (PEO) ₄ propionic acid, succinimidyl ester (3-(azidotetra(ethyleneoxy))propionic acid, succinimidyl ester) 1 mg B10184 biotin azide 1 mg B10185 biotin alkyne 1 mg B10188 iodoacetamide azide 1 mg B10189 iodoacetamide alkyne 5-isomer 488 azide (Oregon Green* 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt) 0.5 mg D10181 Oregon Green* 488 alkyne *6-isomer* 0.5 mg T10182 tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine 5-carboxamido-(6-azidohexanyl)) *5-isomer* 0.5 mg Related Products A10044 EdU (5-ethynyl-2'-deoxyuridine) .50 mg C10102 Click-iT* AHA (L-azidohomoalanine) *for nascent protein synthesis* 5 mg
A10270 Alexa Fluor® 594 azide (Alexa Fluor® 594 carboxamido-(6-azidohexanyl), bis(triethylammonium salt)) 0.5 mg A10275 Alexa Fluor® 594 alkyne (Alexa Fluor® 594 carboxamido-(5-(and 6-)propargyl), bis(triethylammonium salt)) 0.5 mg A10277 Alexa Fluor® 647 azide, triethylammonium salt 0.5 mg A10278 Alexa Fluor® 647 alkyne, triethylammonium salt 0.5 mg A10279 alkyne, succinimidyl ester (3-propargyloxypropanoic acid, succinimidyl ester) 1 mg A10280 azido (PEO)₄ propionic acid, succinimidyl ester (3-(azidotetra(ethyleneoxy))propionic acid, succinimidyl ester) 1 mg B10184 biotin azide 1 mg B10185 biotin alkyne 1 mg B10188 iodoacetamide azide 1 mg B10189 iodoacetamide alkyne 1 mg D10180 Oregon Green® 488 azide (Oregon Green® 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt) 0.5 mg D10181 Oregon Green® 488 alkyne *6-isomer* 0.5 mg T10182 tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine, propargylamide) *5-isomer* 0.5 mg Related Products A10044 EdU (5-ethynyl-2'-deoxyuridine) 50 mg C10102 Click-iT® AHA (L-azidohomoalanine) *for nascent protein synthesis* 5 mg
A10275 Alexa Fluor® 594 alkyne (Alexa Fluor® 594 carboxamido-(5-(and 6-)propargyl), bis(triethylammonium salt)) 0.5 mg A10277 Alexa Fluor® 647 azide, triethylammonium salt 0.5 mg A10278 Alexa Fluor® 647 alkyne, triethylammonium salt 0.5 mg A10279 alkyne, succinimidyl ester (3-propargyloxypropanoic acid, succinimidyl ester) 1 mg A10280 azido (PEO)4 propionic acid, succinimidyl ester (3-(azidotetra(ethyleneoxy))propionic acid, succinimidyl ester) 1 mg B10184 biotin azide 1 mg B10185 biotin alkyne 1 mg I10188 iodoacetamide azide 1 mg I10189 iodoacetamide alkyne 1 mg I10180 Oregon Green® 488 azide (Oregon Green® 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt) 0.5 mg O10181 Oregon Green® 488 alkyne *6-isomer* 0.5 mg T10182 tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine 5-carboxamido-(6-azidohexanyl)) *5-isomer* 0.5 mg T10183 tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer* 0.5 mg Related Products A10044 EdU (5-ethynyl-2'-deoxyuridine) 50 mg C10102 Click-iT® AHA (L-azidohomoalanine) *for nascent protein synthesis* 5 mg
A10277 Alexa Fluor® 647 azide, triethylammonium salt
A10278 Alexa Fluor® 647 alkyne, triethylammonium salt 0.5 mg A10279 alkyne, succinimidyl ester (3-propargyloxypropanoic acid, succinimidyl ester) 1 mg A10280 azido (PEO) ₄ propionic acid, succinimidyl ester (3-(azidotetra(ethyleneoxy)))propionic acid, succinimidyl ester) 1 mg B10184 biotin azide 1 mg B10185 biotin alkyne 1 mg I10188 iodoacetamide azide 1 mg I10189 iodoacetamide alkyne 1 mg O10180 Oregon Green® 488 azide (Oregon Green® 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt) 0.5 mg O10181 Oregon Green® 488 alkyne *6-isomer* 0.5 mg T10182 tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine 5-carboxamido-(6-azidohexanyl)) *5-isomer* 0.5 mg T10183 tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer* 0.5 mg Related Products A10044 EdU (5-ethynyl-2'-deoxyuridine) 50 mg C10102 Click-iT® AHA (L-azidohomoalanine) *for nascent protein synthesis* 5 mg
A10279 alkyne, succinimidyl ester (3-propargyloxypropanoic acid, succinimidyl ester). 1 mg A10280 azido (PEO) ₄ propionic acid, succinimidyl ester (3-(azidotetra(ethyleneoxy))propionic acid, succinimidyl ester) 1 mg B10184 biotin azide. 1 mg B10185 biotin alkyne. 1 mg I10188 iodoacetamide azide. 1 mg I10189 iodoacetamide alkyne. 1 mg O10180 Oregon Green* 488 azide (Oregon Green* 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt) 0.5 mg O10181 Oregon Green* 488 alkyne *6-isomer* 0.5 mg T10182 tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine 5-carboxamido-(6-azidohexanyl)) *5-isomer* 0.5 mg T10183 tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer* 0.5 mg Related Products A10044 EdU (5-ethynyl-2'-deoxyuridine) 50 mg C10102 Click-iT* AHA (L-azidohomoalanine) *for nascent protein synthesis* 5 mg
A10280 azido (PEO) ₄ propionic acid, succinimidyl ester (3-(azidotetra(ethyleneoxy))propionic acid, succinimidyl ester) 1 mg B10184 biotin azide 1 mg B10185 biotin alkyne 1 mg I10188 iodoacetamide azide 1 mg I10189 iodoacetamide alkyne 1 mg O10180 Oregon Green® 488 azide (Oregon Green® 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt) 0.5 mg O10181 Oregon Green® 488 alkyne *6-isomer* 0.5 mg T10182 tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine 5-carboxamido-(6-azidohexanyl)) *5-isomer* 0.5 mg T10183 tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer* 0.5 mg Related Products A10044 EdU (5-ethynyl-2'-deoxyuridine) 50 mg C10102 Click-iT® AHA (L-azidohomoalanine) *for nascent protein synthesis* 5 mg
B10184biotin azide1 mgB10185biotin alkyne1 mgI10188iodoacetamide azide1 mgI10189iodoacetamide alkyne1 mgO10180Oregon Green® 488 azide (Oregon Green® 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt)0.5 mgO10181Oregon Green® 488 alkyne *6-isomer*0.5 mgT10182tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine 5-carboxamido-(6-azidohexanyl)) *5-isomer*0.5 mgT10183tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer*0.5 mgRelated ProductsA10044EdU (5-ethynyl-2'-deoxyuridine)50 mgC10102Click-iT® AHA (L-azidohomoalanine) *for nascent protein synthesis*5 mg
B10185biotin alkyne.1 mgI10188iodoacetamide azide1 mgI10189iodoacetamide alkyne.1 mgO10180Oregon Green* 488 azide (Oregon Green* 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt)0.5 mgO10181Oregon Green* 488 alkyne *6-isomer*0.5 mgT10182tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine 5-carboxamido-(6-azidohexanyl)) *5-isomer*0.5 mgT10183tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer*0.5 mgRelated ProductsA10044EdU (5-ethynyl-2'-deoxyuridine)50 mgC10102Click-iT* AHA (L-azidohomoalanine) *for nascent protein synthesis*5 mg
I10188iodoacetamide azide1 mgI10189iodoacetamide alkyne1 mgO10180Oregon Green® 488 azide (Oregon Green® 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt)0.5 mgO10181Oregon Green® 488 alkyne *6-isomer*0.5 mgT10182tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine 5-carboxamido-(6-azidohexanyl)) *5-isomer*0.5 mgT10183tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer*0.5 mgRelated Products410044EdU (5-ethynyl-2'-deoxyuridine)50 mgC10102Click-iT® AHA (L-azidohomoalanine) *for nascent protein synthesis*5 mg
I10189iodoacetamide alkyne1 mgO10180Oregon Green® 488 azide (Oregon Green® 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt)0.5 mgO10181Oregon Green® 488 alkyne *6-isomer*0.5 mgT10182tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine 5-carboxamido-(6-azidohexanyl)) *5-isomer*0.5 mgT10183tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer*0.5 mgRelated Products8A10044EdU (5-ethynyl-2'-deoxyuridine)50 mgC10102Click-iT® AHA (L-azidohomoalanine) *for nascent protein synthesis*5 mg
O10180 Oregon Green® 488 azide (Oregon Green® 488 6-carboxamido-(6-azidohexanyl), triethylammonium salt). O10181 Oregon Green® 488 alkyne *6-isomer*. T10182 tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine 5-carboxamido-(6-azidohexanyl)) *5-isomer*. 0.5 mg T10183 tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer*. 0.5 mg Related Products A10044 EdU (5-ethynyl-2'-deoxyuridine). 50 mg C10102 Click-iT® AHA (L-azidohomoalanine) *for nascent protein synthesis*. 5 mg
O10181 Oregon Green* 488 alkyne *6-isomer*. 0.5 mg T10182 tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine 5-carboxamido-(6-azidohexanyl)) *5-isomer*. 0.5 mg T10183 tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer* 0.5 mg Related Products A10044 EdU (5-ethynyl-2'-deoxyuridine) 50 mg C10102 Click-iT* AHA (L-azidohomoalanine) *for nascent protein synthesis* 5 mg
T10182tetramethylrhodamine (TAMRA) azide (tetramethylrhodamine 5-carboxamido-(6-azidohexanyl)) *5-isomer*.0.5 mgT10183tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer*.0.5 mgRelated Products8dU (5-ethynyl-2'-deoxyuridine).50 mgC10102Click-iT* AHA (L-azidohomoalanine) *for nascent protein synthesis*.5 mg
T10183tetramethylrhodamine (TAMRA) alkyne (5-carboxytetramethylrhodamine, propargylamide) *5-isomer*0.5 mgRelated ProductsA10044EdU (5-ethynyl-2'-deoxyuridine)50 mgC10102Click-iT* AHA (L-azidohomoalanine) *for nascent protein synthesis*5 mg
Related ProductsA10044EdU (5-ethynyl-2'-deoxyuridine)50 mgC10102Click-iT* AHA (L-azidohomoalanine) *for nascent protein synthesis*5 mg
A10044 EdU (5-ethynyl-2'-deoxyuridine) 50 mg C10102 Click-iT® AHA (L-azidohomoalanine) *for nascent protein synthesis* 5 mg
C10102 Click-iT® AHA (L-azidohomoalanine) *for nascent protein synthesis*
C10186 Click-iT® HPG (L-homopropargylglycine) *for nascent protein synthesis*
C10248 Click-iT® farnesyl alcohol, azide *mixed isomers*
C10249 Click-iT® geranylgeranyl alcohol, azide *mixed isomers*
C10264 Click-iT® fucose alkyne (tetraacetyl fucose alkyne)
C10265 Click-iT® palmitic acid, azide (15-azidopentadecanoic acid)
C10268 Click-iT® myristic acid, azide (12-azidododecanoic acid)
C10269 Click-iT® Cell Reaction Buffer Kit. 1 kit
C10276 Click-iT® Protein Reaction Buffer Kit
C33365 Click-iT® GalNAz metabolic glycoprotein labeling reagent (tetraacetylated <i>N</i> -azidoacetylgalactosamine) *for <i>O</i> -linked glycoproteins*
5.2 mg
C33366 Click-iT® ManNAz metabolic glycoprotein labeling reagent (tetraacetylated <i>N</i> -azidoacetyl- <i>D</i> -mannosamine)
for sialic acid glycoproteins *5.2 mg*
C33367 Click-iT® GlcNAz metabolic glycoprotein labeling reagent (tetraacetylated <i>N</i> -azidoacetylglucosamine)
for O-GlcNAC-modified proteins *5.2 mg*
C33368 Click-iT® O-GlcNAc Enzymatic Labeling System *for O-linked GlcNAc glycoproteins* *10 labelings*
C33370 Click-iT® Tetramethylrhodamine (TAMRA) Protein Analysis Detection Kit *UV/532 nm excitation* *10 reactions*
C33371 Click-iT® Dapoxyl® Protein Analysis Detection Kit *for UV excitation* *10 reactions*
C33372 Click-iT® Biotin Protein Analysis Detection Kit *10 reactions*
E10187 EdU (5-ethynyl-2'-deoxyuridine) 500 mg

Contact Information

Molecular Probes, Inc.

29851 Willow Creek Road Eugene, OR 97402 Phone: (541) 465-8300 Fax: (541) 335-0504

Customer Service:

6:00 am to 4:30 pm (Pacific Time) Phone: (541) 335-0338 Fax: (541) 335-0305 probesorder@invitrogen.com

Toll-Free Ordering for USA:

Order Phone: (800) 438-2209 Order Fax: (800) 438-0228

Technical Service:

8:00 am to 4:00 pm (Pacific Time) Phone: (541) 335-0353 Toll-Free (800) 438-2209 Fax: (541) 335-0238 probestech@invitrogen.com

Invitrogen European Headquarters

Invitrogen, Ltd. 3 Fountain Drive Inchinnan Business Park Paisley PA4 9RF, UK Phone: +44 (0) 141 814 6100 Fax: +44 (0) 141 814 6260 Email: euroinfo@invitrogen.com Technical Services: eurotech@invitrogen.com

For country-specific contact information, visit www.invitrogen.com.

Further information on Molecular Probes products, including product bibliographies, is available from your local distributor or directly and the product of the product bibliographies and the product of the productfrom Molecular Probes. Customers in Europe, Africa and the Middle East should contact our office in Paisley, United Kingdom. All others should contact our Technical Service Department in Eugene, Oregon.

Molecular Probes products are high-quality reagents and materials intended for research purposes only. These products must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Please read the Material Safety Data Sheet provided for each product; other regulatory considerations may apply.

Limited Use Label License No. 223: Labeling and Detection Technology

The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise transfer (b) its components or (c) materials made using this product or its components or (c) materials made using this product or its components or (d) materials made using this product or its components or (d) materials made using the product or its components or (d) materials made using the product or (d) materials made using the product of the party of the product of t erwise use this product or its components or materials made using this product or its components for Commercial Purposes. The buyer may transfer information or materials made through the use of this product to a scientific collaborator, provided that such transfer is not for any Commercial Purpose, and that such collaborator agrees in writing (a) to not transfer such materials to any third party, and (b) to use such transferred materials and/or information solely for research and not for Commercial Purposes. Commercial Purposes means any activity by a party for consideration and may include, but is not limited to: (1) use of the product or its components in manufacturing; (2) use of the product or its components to provide a service, information, or data; (3) use of the product or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product or its components, whether or not such product or its components are resold for use in research. Invitrogen Corporation will not assert a claim against the buyer of infringement of the above patents based upon the manufacture, use or sale of a therapeutic, clinical diagnostic, vaccine or prophylactic product developed in research by the buyer in which this product or its components was employed, provided that neither this product nor any of its components was used in the manufacture of such product. If the purchaser is not willing to accept the limitations of this limited use statement, Invitrogen is willing to accept return of the product with a full refund. For information on purchasing a license to this product for purposes other than research, contact Molecular Probes, Inc., Business Development, 29851 Willow Creek Road, Eugene, OR 97402, Tel: (541) 465-8300. Fax: (541) 335-0354.

Several Molecular Probes products and product applications are covered by U.S. and foreign patents and patents pending. All names containing the designation * are registered with the U.S. Patent and Trademark Office.

Copyright 2009, Molecular Probes, Inc. All rights reserved. This information is subject to change without notice.