

Application Guide

IRDye[®] Peptide Labeling



Published April 2013. Revised April 2022. The most recent version of this document is posted at [licor.com/bio/support](https://www.licor.com/bio/support).

Table of Contents

	Page
I. Introduction	3
II. Peptide Labeling Options	4
III. Labeling Reaction Techniques	5
Aqueous Solution-Phase Labeling	5
Organic Solution-Phase Labeling	6
Purification	7
Quantification	7
IV. Examples	9
V. References	9

I. Introduction

Peptides labeled with fluorescent dyes are important as probes for *in vivo* imaging and as substrates for enzyme activity assays. Near-infrared (NIR) fluorophores such as IRDye® 800CW (excitation_{Max} 774 nm; emission_{Max} 789 nm) can offer improved sensitivity because of low NIR autofluorescence from tissues, cells, biological materials, or drug compounds.

Fluorophores can also be combined with appropriate quencher dyes to create fluorogenic peptide probes, such as those used for protease activity detection (1 – 4). IRDye QC-1 can efficiently quench a wide range of fluorophores spanning the visible and NIR spectra (~500–800 nm) in a fluorescence resonance energy transfer (FRET) system. Together, IRDye 800CW and IRDye QC-1 comprise an optimal fluorophore-quencher pair for incorporation into NIR fluorogenic peptides (Figure 1).

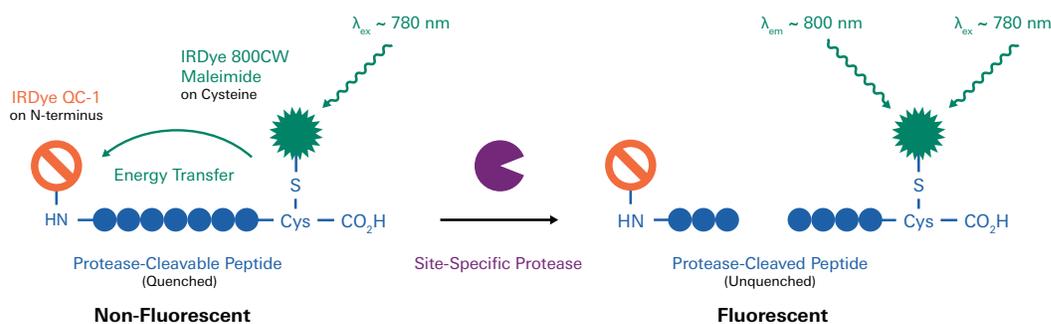


Figure 1. NIR fluorogenic peptide probe using IRDye 800CW donor and IRDye QC-1 quencher.

IRDye fluorophores and quenchers are amenable to a wide variety of conjugation reactions that yield stable, covalently-labeled peptides. Furthermore, site-specific dye labeling can be achieved by employing orthogonally reactive groups in peptide design. Because IRDye labels have been optimized for detection in aqueous biological environments, conjugation reactions are best performed in aqueous-phase.

II. Peptide Labeling Options

IRDye NHS esters react with unprotected amines in peptides. NHS ester reactions usually proceed quickly and cleanly in either organic or aqueous solvents. While NHS esters continue to be widely used tools for biomolecule modification, their application to large peptides may be complicated by factors, such as steric hindrance and multiple labeling sites.

To address the limitations of NHS ester chemistry, orthogonal technologies such as IRDye Maleimides and Click Chemistry Reagents have been developed to label non-amine functional groups. Orthogonal chemistry can also facilitate dual- or multi-dye labeling. As with NHS ester chemistry, optimal non-amine labeling is usually achieved by aqueous solution-phase reactions. Table 1 summarizes general strategies for labeling peptides with IRDye reagents.

Table 1. Appropriate IRDye reagents for labeling peptides

Desired Peptide Labeling Site	IRDye Reagent	Notes
Unblocked N-terminus or Side-Chain Amine (e.g., lysine)	NHS Ester	Perform in mildly basic aqueous solutions; peptides containing >1 amine may yield product mixtures
Sulfhydryl (e.g., cysteine)	Maleimide	Perform in mildly acidic aqueous solutions for selective labeling in the presence of amines
Alkyne	Azide	Copper-catalyzed Click Chemistry for selective labeling in the presence of amines
Azide	Alkyne	Copper-catalyzed Click Chemistry for selective labeling in the presence of amines
Azide	DBCO	Copper-free Click Chemistry for selective labeling in the presence of amines
DBCO	Azide	Copper-free Click Chemistry for selective labeling in the presence of amines

IRDye Maleimides are designed to react selectively with sulfhydryl groups in the presence of amines. Optimal selectivity is usually obtained with labeling reactions performed in mildly acidic (pH 6.5) aqueous solvents. At acidic pH, amine reactivity is greatly reduced because they are prevalently protonated. Sulfhydryl groups can be installed in peptides with cysteine or commercially available thiol-linkers.

Note: Peptides containing sulfhydryls can spontaneously form inactive dimers or cyclize internally. Therefore, it may be necessary to pre-treat the peptide with a reducing agent such as tris(2-carboxyethyl)- phosphine (TCEP) to reveal reactive sulfhydryl functional groups.

IRDye Click Chemistry reagents are mostly inert to naturally occurring functional groups such as amines and hydroxyls. Click Chemistry utilizes pairs of functional groups that can react either in the presence (5) or absence (6) of a copper catalyst. Click Chemistry functional groups can also be readily installed in peptides with commercially available reagents.

III. Labeling Reaction Techniques

Aqueous Solution-Phase Labeling

For all labeling reactions, it is critical to omit components that can interfere with the coupling of the peptide and the dye. Peptides should be dissolved in aqueous buffers that do not contain extraneous nucleophiles such as Tris, sodium azide, DTT, BME, etc.

To maximize the labeling efficiency, the peptide should be dissolved at the highest practical concentration that still maintains a homogeneous reaction mixture. The aqueous reaction mixture can be directly purified by reverse-phase HPLC (see "Purification" on page 7). Table 2 provides general parameters for aqueous solution-phase labeling of peptides with IRDye reagents.

Table 2. General parameters for labeling peptides in aqueous solutions

IRDye Reagent	Dye Equivalents per Labeling Site	Reaction Buffer (pH)	Reaction Temperature	Reaction Time (hours)
NHS Ester	1–2	Phosphate (7.5–8.5)	4 °C to 37 °C	1–3 for ambient to 37 °C or overnight for 4 °C
Maleimide	1–2	Phosphate (6.5–7.5)	4 °C to 37 °C	2–3
Alkynyl†	1	Variable (6.5 to 8.5)	Ambient to 37 °C	1–4
Azide†,*	1	Variable (6.5 to 8.5)	Ambient to 37 °C	1–4
DBCO*	1	Variable (6.5 to 8.5)	Ambient to 37 °C	1–4

* Copper-catalyzed Click Chemistry (5)

† Copper-free Click Chemistry (6)

Organic Solution-Phase Labeling

Certain peptides, such as those comprised mainly of hydrophobic amino acids, may not be amenable to aqueous solution-phase labeling. In these atypical cases, IRDye reagents can be used for organic solution-phase labeling; however, this reaction may be slower than its aqueous counterpart.

Again, peptides should be dissolved in solvents devoid of extraneous nucleophiles such as Tris, sodium azide, DTT, BME, etc. The optimal reaction solvent is anhydrous dimethyl sulfoxide (DMSO). When labeling with IRDye NHS esters and Maleimides, including a tertiary amine base such as *N,N*-diisopropylethylamine (DIPEA) is necessary to promote the reaction. After the reaction has completed, the crude IRDye labeled peptide can be precipitated from the DMSO solution by dropwise addition into anhydrous diethyl ether.

IRDye labels impart hydrophilicity and can facilitate the purification of hydrophobic peptides by reverse-phase HPLC (see "Purification" on the facing page). Table 3 provides general parameters for organic solution-phase labeling of peptides with IRDye reagents.

Table 3. General parameters for labeling peptides in organic solutions

IRDye Reagent	Dye Equivalents per Labeling Site	Base Equivalents per Labeling Site	Reaction Temperature	Reaction Time (hours)
NHS Ester	1–2	2–4	Ambient to 37 °C	2–12
Maleimide	1–2	0–2	Ambient to 37 °C	2–12
Alkyne†	1	None	Ambient to 37 °C	1–4
Azide*	1	None	Ambient to 37 °C	1–4
DBCO*	1	None	Ambient to 37 °C	1–4

* Copper-Catalyzed Click Chemistry ([5](#))

† Copper-Free Click Chemistry ([6](#))

Purification

Reverse-phase HPLC purification delivers IRDye labeled peptides with the highest purity. The recommended mobile phase is a gradient mixture of acetonitrile and water buffered with triethylammonium acetate (TEAA, 50 mM, pH 6.0). The TEAA provides ion-pairing for better retention behavior of the hydrophilic IRDye labeled peptides on the column.

As a consequence, the predominant counter-ion of the purified peptide will be triethylammonium, which may interfere with downstream biological experiments. Prior to lyophilization, the purified peptide should be ion-exchanged by eluting through an appropriate resin (e.g., Amberlite™) or by dialysis.

For flexibility in development and/or troubleshooting, the HPLC system should be equipped with a diode array detector (DAD) and be able to monitor at the absorption maxima of all dyes used.

Quantification

UV-Vis absorbance spectroscopy is the best method for quantifying the amount of purified IRDye labeled peptide. To determine the concentration of a stock solution containing a singly-labeled peptide, dilute an aliquot of the stock solution in 1X PBS, measure the absorbance spectra at the dye-specific maximum and use the following equation:

$$\text{Conc of IRDye peptide (mg/mL)} = \frac{A_{\text{IRDye}}}{\epsilon_{\text{IRDye}}} \times \text{Dilution Factor} \times \text{MW}_{\text{IRDye peptide}}$$

In which:

- A_{IRDye} is the measured absorbance at the dye-specific maximum (See Table 4)
- ϵ_{IRDye} is the extinction coefficient for the dye in 1X PBS (See Table 4)
- Dilution Factor is the fold dilution of the IRDye peptide in 1X PBS
- $\text{MW}_{\text{IRDye peptide}}$ is the molecular weight of the IRDye peptide

Table 4. Parameters for UV-Vis quantification of IRDye labeled peptides in 1X PBS

Label	Dye Maximum (nm) for Measuring Absorbance	ε _{IRDye}
IRDye 800CW	778	240,000
IRDye 800RS	767	200,000
IRDye 750	756	260,000
IRDye 680RD	672	165,000
IRDye 680LT	676	250,000
IRDye 650	648	230,000
IRDye QC-1	737	96,000

For a pure stock solution of a FRET-quenched peptide labeled with exactly one IRDye 800CW and one IRDye QC-1: Determine the concentration of a stock solution by diluting an aliquot of the stock solution in methanol, measuring the absorbance spectra at 778 nm and 850 nm and using the following equation:

$$\text{Conc of IRDye 800CW QC-1 peptide (mg/mL)} = \frac{A_{778} - (1.265 \times A_{850})}{300,000} \times \text{Dilution Factor} \times \text{MW}_{\text{IRDye 800CW QC-1 peptide}}$$

In which:

- A_{778} is the **measured absorbance** at the maximum for IRDye 800CW in methanol
- 300,000 is the extinction coefficient for IRDye 800CW in methanol
- 1.265 is the correction factor for the ratio of A_{778}/A_{850} for IRDye QC-1 in methanol
- A_{850} is the **measured absorbance** at the maximum for IRDye QC-1 in methanol
- Dilution Factor is the fold dilution of the IRDye 800CW QC-1-peptide in methanol
- $\text{MW}_{\text{IRDye 800CW QC-1 peptide}}$ is the molecular weight of the IRDye 800CW QC-1 peptide

IV. Examples

- **For an example of aqueous solution-phase labeling with IRDye 800CW NHS ester, see** Davies-Venn, C.A., Angermiller, B., Wilganowski, N., Ghosh, P., Harvey, B.R., Wu, G., et al. (2012). Albumin-Binding Domain Conjugate for Near-Infrared Fluorescence Lymphatic Imaging. *Mol. Imaging Biol.*, 14(3), 301–14. [DOI: 10.1007/s11307-011-0499-x](https://doi.org/10.1007/s11307-011-0499-x)
- **For an example of organic solution-phase labeling with IRDye 800CW NHS ester, see** Chen, Y., Dhara, S., Banerjee, S.R., Byun, Y., Pullambhatla, M., Mease, R.C., and Pomper, M. (2009). A low molecular weight PSMA-based fluorescent imaging agent for cancer. *Biochem. Biophys. Res. Commun.*, 390(3), 624–9. [DOI: 10.1016/j.bbrc.2009.10.017](https://doi.org/10.1016/j.bbrc.2009.10.017)
- **For an example of aqueous solution-phase labeling with IRDye 800CW Maleimide, see** Ye, Y., Zhu, L., Ma, Y., Niu, G., and Chen, X. (2011). Synthesis and evaluation of new iRGD peptide analogs for tumor optical imaging. *Bioorg. Med. Chem. Lett.*, 21(4), 1146–50. [DOI: 10.1016/j.bmcl.2010.12.112](https://doi.org/10.1016/j.bmcl.2010.12.112)
- **For examples of dual-labeling with IRDye QC-1 NHS ester and various fluorophores, see** Sun, X., Zhang, A., Baker, B., Sun, L., Howard, A., Buswell, J., et al. (2011). Development of SNAP-Tag Fluorogenic Probes for Wash-Free Fluorescence Imaging. *ChemBioChem*, 12(14), 2217–26. [DOI: 10.1002/cbic.201100173](https://doi.org/10.1002/cbic.201100173)

V. References

1. Matayoshi, E.D., Wang, G.T., Krafft, G.A., and Erickson, J. (1990). Novel fluorogenic substrates for assaying retroviral proteases by resonance energy transfer. *Science*, 247(4945), 954–8. [DOI: 10.1126/science.2106161](https://doi.org/10.1126/science.2106161)
2. Bullok, K.E., Maxwell, D., Kesarwala, A.H., Gammon, S., Prior, J.L., Snow, M., et al. (2007). Biochemical and *in vivo* characterization of a small, membrane-permeant, caspase-activatable far-red fluorescent peptide for imaging apoptosis. *Biochem.*, 46(13), 4055–65. [DOI: 10.1021/bi061959n](https://doi.org/10.1021/bi061959n)
3. Zheng, G., Chen, J., Stefflova, K., Jarvi, M., Li, H., and Wilson, B.C. (2007). Photodynamic molecular beacon as an activatable photosensitizer based on protease controlled singlet oxygen quenching and activation. *Proc. Natl. Acad. Sci. USA*, 104 (21), 8989–94. [DOI: 10.1073/pnas.0611142104](https://doi.org/10.1073/pnas.0611142104)
4. Blum, G., Mullins, S.R., Keren, K., Fonovic, M., Jedeszko, C., Rice, M.J., et al. (2005). Dynamic imaging of protease activity with fluorescently quenched activity-based probes. *Nature Chem. Biol.*, 1(4), 203–9. [DOI: 10.1038/nchembio728](https://doi.org/10.1038/nchembio728)
5. Best, M.D. (2009). Click Chemistry and bioorthogonal reactions: Unprecedented selectivity in the labeling of biological molecules. *Biochemistry*, 48(28), 6571–84. [DOI: 10.1021/bi9007726](https://doi.org/10.1021/bi9007726)
6. Simon, M., Zangemeister-Wittke, U., and Plückthun, A. (2012). Facile double-functionalization of designed ankyrin repeat proteins using Click and thiol chemistries. *Bioconjugate Chem.*, 23(2), 279–86. [DOI:10.1021/bc200591x](https://doi.org/10.1021/bc200591x)



© 2022 LI-COR, Inc. LI-COR and IRDye are trademarks or registered trademarks of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners. Use of IRDye infrared dyes in copper-catalyzed click chemistry reactions is covered by one or more U.S. or foreign patents owned by Scripps Research Institute. For patent information, visit licor.com/patents.

LI-COR Biosciences

4647 Superior Street
Lincoln, NE 68504
Phone: +1-402-467-0700
Toll free: 800-645-4267
biosales@licor.com
licor.com/bio

Regional Offices

LI-COR Biosciences GmbH
Siemensstraße 25A
61352 Bad Homburg
Germany
Phone: +49 (0) 6172 17 17 771
bio-eu@licor.com

LI-COR Biosciences UK Ltd.
St. John's Innovation Centre
Cowley Road • Cambridge
CB4 0WS • United Kingdom
Phone: +44 (0) 1223 422104
bio-eu@licor.com