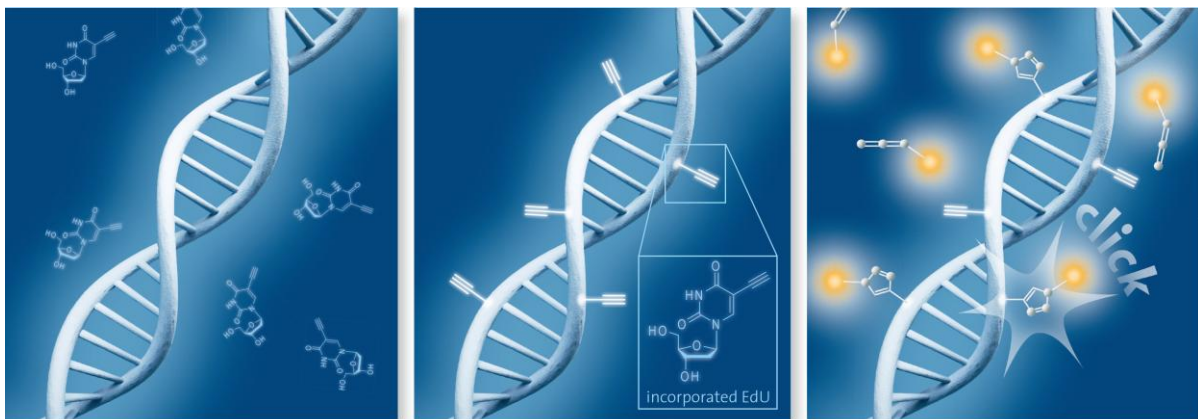




## EdU Flow Cytometry Kit

### User Manual



**Ordering information:***(for detailed kit content see Table 2)***EdU Flow Cytometry Kits for 50 assays:**

Product number	EdU	Used fluorescent dye
BCK-FC488-50	10 mg	6-FAM Azide
BCK-FC555-50	10 mg	5-TAMRA-PEG3-Azide
BCK-FC594-50	10 mg	5/6-Sulforhodamine 101-PEG3-Azide
BCK-FC647-50	10 mg	Cyanine 5 Azide

**EdU Flow Cytometry Kits for 100 assays:**

Product number	EdU	Used fluorescent dye
BCK-FC488-100	20 mg	6-FAM Azide
BCK-FC555-100	20 mg	5-TAMRA-PEG3-Azide
BCK-FC594-100	20 mg	5/6-Sulforhodamine 101-PEG3-Azide
BCK-FC647-100	20 mg	Cyanine 5 Azide

To place your order, please contact us under:

- phone: +49 8158 903867
- fax: +49 8158 903894
- email: [orders@baseclick.eu](mailto:orders@baseclick.eu)

---

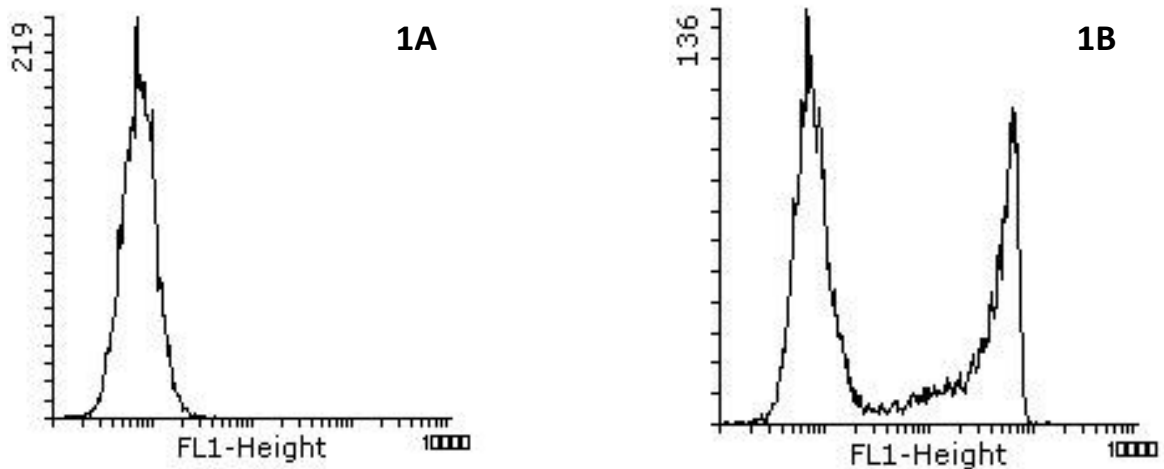
## EdU Flow Cytometry Kit

### Introduction and product description:

The detection of cell proliferation is of utmost importance for assessing cell health, determining genotoxicity or evaluating anticancer drugs. This is normally performed by adding nucleoside analogs like [<sup>3</sup>H]thymidine or 5-bromo-2'-deoxyuridine (BrdU) to cells during replication, and their incorporation into DNA is detected or visualized by autoradiography or with an anti-BrdU-antibody respectively. Both methods exhibit several limitations. Working with [<sup>3</sup>H]thymidine is troublesome because of its radioactivity. Autoradiography is slow and thus not suitable for rapid high-throughput studies. The major disadvantage of BrdU staining is that the double-stranded DNA blocks the access of the anti-BrdU antibody to BrdU units. Therefore samples have to be subjected to harsh denaturing conditions resulting in degradation of the structure of the specimen.

The baseclick *EdU Flow Cytometry Kits* overcome these limitations, providing a superior alternative to BrdU and [<sup>3</sup>H]thymidine assays for directly measuring DNA synthesis. EdU (5-ethynyl-2'-deoxyuridine) is a nucleoside analog to thymidine and is incorporated into DNA during active DNA synthesis. In contrast to BrdU assays, the *EdU Flow Cytometry Assays* are not antibody based and therefore do not require DNA denaturation for detection of the incorporated nucleoside. Instead, the *EdU Flow Cytometry Kits* utilize click chemistry for detection in a variety of dye fluorescent readouts. Furthermore, the streamlined detection protocol reduces both the total number of steps and significantly decreases the total amount of time. The simple click chemistry detection procedure is complete within 30 minutes and is compatible with multiplexing for content and context-rich results.

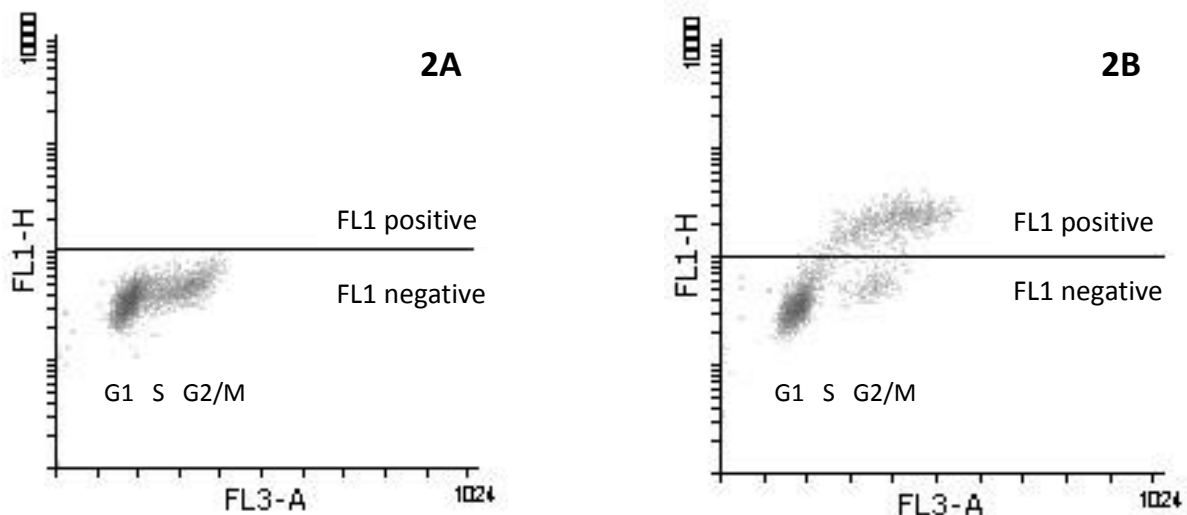
Standard flow cytometry methods are used to determine the percentage of S-phase cells in the population (**Figure 1**).



**Figure 1: Fluorescence histograms of EdU-incorporation with baseclick EdU Flow Cytometry Kit.**

Samples of HeLa cells treated without (**1A**) or with EdU (**1B**) were incubated with 10  $\mu$ M EdU for 2 hours. The click reaction using 6-FAM Azide was performed according to the recommended staining protocol. Fluorescence intensity of 10.000 cells was measured by flow cytometry. The results are presented in form of histograms, showing the cell number in the y-axis and the FL1-Fluorescence in the x-axis. FL1 voltage setting was adjusted according to the fluorescence signal of the negative cell population (333 V with 6-FAM). **1A** represents the negative control of proliferating and non-proliferating cells without EdU incorporation. **1B** shows non-proliferating cells without EdU incorporation (left peak) and proliferating cells (S phase) which have incorporated EdU and are labelled with 6-FAM Azide (right peak).

The baseclick *EdU Flow Cytometry Kit* is compatible with several cell cycle dyes. An example using 6-FAM Azide is illustrated in **Figure 2**.



**Figure 2: Density blots of Propidium Iodide (PI) stained samples.**

Samples of HeLa cells treated without (**2A**) or with EdU (**2B**) were incubated with 10  $\mu$ M EdU for 2 hours. The reaction cocktail carrying 6-FAM Azide was used. After the click reaction DNA was stained using PI (FL3 fluorescent channel). The y-axis presents the FL1-Fluorescence intensity, and the x-axis the content of DNA measured with FL3-area. Cell cycle phases are indicated as G1, S and G2/M phase.

The baseclick *EdU Flow Cytometry Kit* can be used with antibodies against surface and intracellular markers. To ensure the compatibility of your reagent or antibody, please refer to **Table 1**.

**Table 1:** EdU detection dye compatibility

Fluorescent molecule	Compatibility
Organic dyes such as Fluorescein and Alexa dyes	Compatible
PerCP, Allophycocyanin (APC) and APC-based tandems	Compatible
R-phycoerythrin (R-PE) and R-PE based tandems	Use R-PE and R-PE based tandems <b>after</b> the EdU detection reaction
Quantum Dots	Use Quantum Dots <b>after</b> the EdU detection reaction
Fluorescent proteins (e.g. GFP)	Use anti-GFP antibodies* <b>before</b> the EdU detection reaction or use organic dye-based reagents for protein expression detection

\* Compatibility indicates which involved components are unstable in the presence of copper catalyst for the EdU detection reaction (either the fluorescent dye itself or the detection method). Not all GFP antibodies recognize the same antigen site. Rabbit and chicken anti-GFP antibodies result in a good fluorescent amount. The mouse monoclonal antibodies tested are not recommended for this application because they do not generate an acceptable amount of fluorescence.

#### For research use only.

Information in this document is subject to change without notice. baseclick GmbH assumes no responsibility for any errors that may appear in this document.

baseclick GmbH disclaims all warranties with respect to this document, expressed or implied, including but not limited to those of merchantability or fitness for a particular purpose. In no event shall baseclick GmbH be liable, whether in contract, tort, warranty, or under any statute or on any other basis for special, incidental, indirect, punitive, multiple or consequential damages in connection with or arising from this document, including but not limited to the use thereof.

Please read the material safety data sheets (MSDS) provided for each product/component.

#### Cautions:

**DMSO (Component C):** is known to facilitate the entry of organic molecules into tissues. DMSO is irritant and flammable. Handle reagents containing DMSO using equipment and practices appropriate for the hazards posed by such materials. Dispose of the reagents in compliance with all related local arrangements.

**Fixative solution (Component D):** contains paraformaldehyde, which is harmful. Use with appropriate precautions.

**Saponin based permeabilization and wash reagent (Component E):** contains sodium azide, which is highly toxic and yields the extremely toxic hydrazoic acid under acidic conditions. Dilute azide compounds in running water before discarding to avoid accumulation of potentially explosive deposits in plumbing. This solution is orange.

**Literature Citation:** When describing a procedure for publication using this product, please refer to it as *baseclick EdU Flow Cytometry Kit*.

## 1. Materials provided with the Kit and storage conditions

**Table 2:** Contents of the kit and storage conditions

Vial-label	Amount for 50 assays	Amount for 100 assays	Component	Component long term storage	Kit storage
<b>Component A</b>	10 mg	20 mg	5-Ethynyl-deoxyuridine (5-EdU)	-20°C	2 - 8°C Dark Do not freeze Dry
<b>Component B red</b>	130 µL	2 x 130 µL	6-FAM Azide (BCK-FC488) 5-TAMRA-PEG3-Azide (BCK-FC555) 5/6-Sulforhodamine101-PEG3-Azide (BCK-FC594) Cyanine 5 Azide (BCK-FC647)	-20°C dark	
<b>Component C</b>	5 mL	8.5 mL	DMSO	RT	
<b>Component D</b>	5 mL	2 x 5 mL	Fixative solution (4% Paraformaldehyde in PBS)	2 – 8°C	
<b>Component E</b>	50 mL	2 x 50 mL	Saponin-based permeabilization and wash reagent (10x solution)	2 – 8°C	
<b>Component F green</b>	2 mL	2 mL	Catalyst solution	RT	
<b>Component G</b>	400 mg	400 mg	Buffer additive	-20°C	

This kit is stable up to 1 year after receipt, when stored as directed.

## 2. Required Material and Equipment not included in this Kit

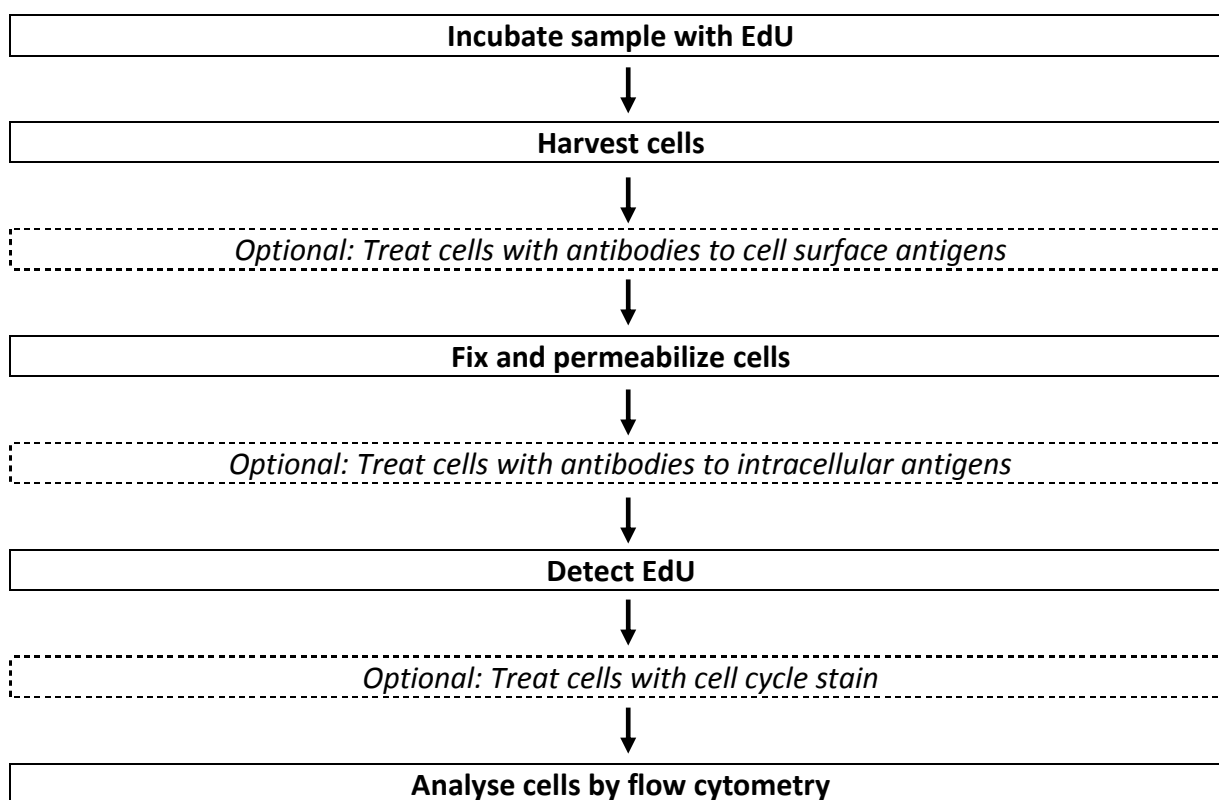
- Adherent cells
- Reaction tubes (size depends on the volume of reaction cocktail needed)
- Buffered saline solution, such as PBS, D-PBS or TBS
- Appropriate cell culture medium
- 1% BSA (bovine serum albumin) in PBS, pH 7.1 – 7.4
- Deionized water or 18 MΩ purified water
- Flow cytometry tubes

## 3. Workflow

The following protocol was developed using an EdU concentration of 10 µM and can be adapted for any cell type. There are many factors which can influence the labeling such as the growth medium, the density and the type of cells. To determine the optimal concentration for your experiment, a range of EdU concentrations should be tested for your cell type and experimental conditions.

Principally, a similar concentration to BrdU can be used for EdU as a starting point. Heparin can be used as anticoagulant for collection, if a whole blood sample is used.

#### Workflow scheme for the *EdU Flow Cytometry Assay*



## 4. Preparation of the stock solutions

- 4.1** Allow all vials to warm to room temperature before opening.
- 4.2** For the preparation of a 10 mM stock solution of EdU, add the appropriate amount of DMSO (**component C**) or aqueous solution (PBS) to EdU (**component A**) according to **table 3** and mix until the compound is completely dissolved. After use, store any remaining solution at -20°C. When stored as directed, this stock solution is stable for up to one year.

**Table 3:** Amounts of DMSO or aqueous solution needed to dissolve EdU to a final concentration of 10 mM

EdU Flow Cytometry Kit	EdU amount	DMSO/aqueous solution amount
50 assays kit	10 mg	4 mL
100 assays kit	20 mg	8 mL

- 4.3** For the preparation of a 10x stock solution of the buffer additive, add 4 mL of deionized water to the **component G** and mix until the compound is dissolved completely. After use, store any remaining solution at -20°C. When stored as directed, this stock solution is stable for up to 6 months. If the solution starts to

develop a brown colour, it has degraded and should be discarded. We recommend preparing aliquots to avoid repeated thaw and freeze cycles!

- 4.4** For the preparation of 500 mL of the 1x saponin-based permeabilization buffer and wash reagent (for 50 assays), add 50 mL of **component E** to 450 mL of 1% BSA in PBS. For the preparation of 1 L of the 1x saponin-based permeabilization buffer and wash reagent (for 100 assays), add 100 mL of **component E** to 900 mL of 1% BSA in PBS. After use, store any remaining solution at 2 - 8°C.

**Note:** The saponin-based permeabilization and wash reagent contains sodium azide. Please see the cautions on page 4.

## 5. Labeling of cells with EdU

- 5.1** Suspend the cells in an appropriate tissue culture medium to obtain optimal cell growth conditions. Please note that the growth of the cells during incubation decelerates, if the temperature changes or the cells are washed prior to incubation with EdU.
- 5.2** For the desired final concentration, add the appropriate amount of EdU to the culture medium and mix well. We recommend using a concentration of 10 µM for 1-2 hours as a starting point. Use higher EdU concentrations for a shorter incubation time. A longer incubation time requires lower EdU concentrations.
- 5.3** The incubation of the cells with EdU should be performed under the optimal conditions for your cell type and for the desired length of time. Various DNA synthesis and proliferation parameters can be evaluated by altering the EdU incubation time or by subjecting the cells to pulse labeling with EdU. Effective time intervals for pulse labeling and the length of each pulse depend on the cell growth rate.
- 5.4** Harvest cells. If performing antibody surface labeling, proceed immediately to step **6**, otherwise continue to step **7**.

## 6. Staining cell-surface antigens with antibodies (optional)

- 6.1** Wash cells with 3 mL of 1% BSA in PBS. Centrifuge to pellet cells and remove supernatant.
- 6.2** Dislodge the pellet and resuspend cells in 1% BSA in PBS at  $1 \times 10^7$  cells/mL.
- 6.3** Add 100 µL of cell suspension or whole blood sample to flow tubes.
- 6.4** Add surface antibodies and mix well.  
**Note:** PE, PE-tandem or Quantum Dot antibody conjugates should not be used before performing the click reaction (step **8**).
- 6.5** Incubate the cells for the recommended length of time and temperature. Protect from light!
- 6.6** Proceed to step **7**.



## 7. Cell fixation and permeabilization

This protocol was developed with a fixation step using 4% Paraformaldehyde in PBS, followed by a saponin-based permeabilization step. The saponin-based permeabilization and wash reagent can be used with cell suspensions containing red blood cells or whole blood as well as with cell suspensions containing different cell types. The morphological light scatter characteristics of leukocytes are maintained by the permeabilization reagent while red blood cells are lysed.

- 7.1** Remove the incubation media and wash the cells with 3 mL of 1% BSA in PBS. Pellet the cells and remove the supernatant.
- 7.2** Dislodge the cell pellet. Add 100  $\mu$ L of the fixative solution (**component D**) to the cells. Mix well and incubate for 15 minutes at room temperature. Protect from light.
- 7.3** Remove the fixation solution and wash the cells with 3 mL of 1% BSA in PBS. Pellet the cells and remove the supernatant. If red blood cells or haemoglobin are present in the sample, repeat the washing step. Remove all residual blood cell debris and haemoglobin before proceeding.
- 7.4** Dislodge the cell pellet. Resuspend the cells in 100  $\mu$ L of 1x saponin-based permeabilization buffer in PBS (prepared in **4.4**). Mix well and proceed to step **8**. for the click reaction.

## 8. EdU detection

- 8.1** Prepare the assay cocktail in the same order as described in **table 4**. If the ingredients are not added in the order listed, the reaction will not proceed optimally or might even fail.

**Important:** Once the assay cocktail is prepared, use it immediately, at least within the next 15 minutes!

**Table 4:** Click assay cocktails

Material	Component	Number of assays				
		1	2	3	5	10
PBS, D-PBS or TBS	Not provided!	438 $\mu$ L	875 $\mu$ L	1.32 mL	2.19 mL	4.38 mL
Catalyst solution	<b>F - green</b>	10 $\mu$ L	20 $\mu$ L	30 $\mu$ L	50 $\mu$ L	100 $\mu$ L
Dye Azide (10 mM)	<b>B - red</b>	2.5 $\mu$ L	5 $\mu$ L	7.5 $\mu$ L	12.5 $\mu$ L	25 $\mu$ L
Buffer additive (10x) (prepared in <b>4.3</b> )	<b>G</b>	50 $\mu$ L	100 $\mu$ L	150 $\mu$ L	250 $\mu$ L	500 $\mu$ L
<b>Total Volume</b>	-	<b>500 <math>\mu</math>L</b>	<b>1 mL</b>	<b>1.5 mL</b>	<b>2.5 mL</b>	<b>5 mL</b>

- 8.2** Add the appropriate amount of the assay cocktail to the cells and mix well to distribute the assay solution evenly.
- 8.3** Incubate the assay mixture for 30 minutes at room temperature. Protect from light!

- 8.4** Wash the cells with 3 mL of 1x saponin based permeabilization and wash reagent (prepared in **4.4**). Pellet the cells and remove the supernatant. Dislodge the cell pellet. If proceeding with intracellular antibody labeling in step **9**, resuspend the cells in 100  $\mu$ L of 1x saponin-based permeabilization and wash reagent. Otherwise, add 500  $\mu$ L of 1x saponin-based permeabilization and wash reagent and proceed with step **10** for analyzing the cells with a flow cytometer.

**Important:** Keep the samples protected from light during the whole procedure.

### **9. Staining intracellular or surface antigens (optional)**

- 9.1** Add antibodies against intracellular antigens or against surface antigens that use RPE, PR-tandem or Quantum Dot antibody conjugates. Mix well.
- 9.2** Incubate the tubes for the time and temperature required for antibody staining. Protect from light.
- 9.3** Wash each tube with 3 mL 1x saponin-based permeabilization and wash reagent (prepared in **4.4**). Pellet the cells and remove the supernatant. Dislodge the cell pellet and resuspend the cells in 500  $\mu$ L of 1x saponin-based permeabilization and wash reagent.
- 9.4** Proceed with step **10** for analyzing the cells with a flow cytometer.

### **10. Imaging and analysis**

Use a low flow rate during acquisition, if a traditional flow cytometer with a hydrodynamic focusing is used to measure the total DNA content. The same collection rate and cell concentration should be used for each sample within an experiment. Detect the fluorescent signal generated by DNA content stains with linear amplification. The fluorescent signal generated by EdU labeling is best detected with logarithmic amplification.

The Excitation and emission maxima of the available dyes are listed in **table 5**.

**Table 5:** Emission and excitation maxima of the available dyes.

Product number	Dye	Excitation (nm)	Emission (nm)	Filter
BCK-FC488	6-FAM Azide	496	516	Green
BCK-FC555	5-TAMRA-PEG3-Azide	546	579	Violet
BCK-FC594	5/6-Sulforhodamine 101-PEG3-Azide	584	603	Orange
BCK-FC647	Cyanine 5 Azide	646	662	Red





baseclick GmbH

Bahnhofstraße 9-15

82327 Tutzing, Germany

Phone: +49 8158 903867

Fax: +49 8158 903894

Email: [info@baseclick.eu](mailto:info@baseclick.eu)