

ReadiUse™ Preactivated PE-iFluor 647 Tandem

Ordering Information

Product Number: 2577 (1mg)

Storage Conditions

Store desiccated at -20 °C and protected from light.
Expiration date is 12 months from the date of receipt.

Introduction:

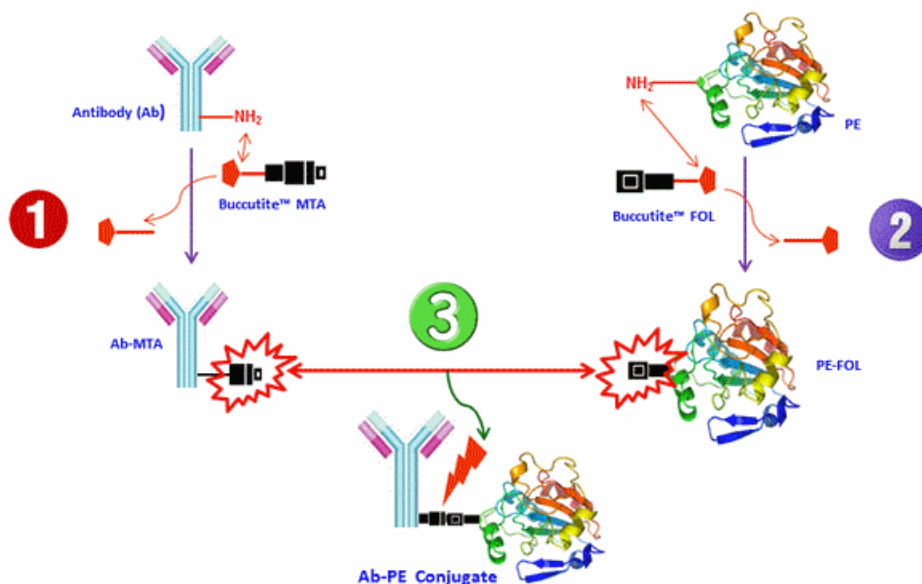
R-Phycoerythrin (PE) is isolated from red algae. Its primary absorption peak is at 565 nm with secondary peaks at 496 and 545 nm. ReadUse™ Preactivated PE-iFluor647 Tandem has been pre-activated with our proprietary linker Buccutite™ FOL, which readily reacts with Buccutite™ MTA-containing molecules under extremely mild neutral conditions. Our Buccutite™ bioconjugation system is much more robust and easier to use and enables faster and quantitative conjugation of biomolecules with higher efficiencies and yields.

Storage Conditions

Store at -20 °C, desiccated and protected from light. Expiration date is one year from the date of receipt.

Conjugation Principle

PE was premodified with our Buccutite™ FOL as shown in Step 2 (provided). Your antibody (or other proteins) is modified with our Buccutite™ MTA to give MTA-modified protein as shown in Step 1. The MTA-modified protein readily reacts with FOL-modified PE (provided) to give the desired PE-antibody conjugate as shown in Step 3.



Conjugation Protocols:

- Pre-activate Antibody with MTA
 - Reconstitute MTA in DMSO at ~10 mg/mL.
Note: Please store unused MTA at -20°C and could be used up to two freeze and thaw cycles.
 - Prepare target antibody (Ab) in pH=8.5~9.0 buffer at concentration above 1mg/ml.
 - Add MTA to Ab solution at the ratio of 8~10ug MTA /100 ug Ab.
 - Mix well and react at RT for 60 minutes, rotating during the reaction.
 - Purify the reaction mixture with desalting column to remove unreacted MTA and exchange buffer to PBS or buffer of your choice.

- 1.6 Collect the MTA-activated Ab, and estimate the concentration by 70% yield of the original starting amount.
2. Conjugate with Pre-activated PE-iFluor647 Tandem
 - 2.1 Reconstitute pre-activated PE- iFluor647 Tandem in 100 µL ddH₂O to 10mg/mL.
Note: Reconstituted pre-activated PE- iFluor647 Tandem is not stable and could be stored at 4°C for one month, please kept from light.
 - 2.2 Add PE-iFluor647 directly to MTA-activated target Ab solution (from 1.5) at the ratio of 300µg PE-iFluor647 /100µg MTA-activated Ab.
 - 2.3 Rotate the mixture for 1~2 hours at room temperature
 - 2.4 The Ab/ PE-iFluor647 conjugate is now ready to use.
Optional: Ab/ PE-iFluor647 Conjugate could be further purified through size exclusion chromatography to get best performance.
Note1: The antibody conjugate should be stored at > 0.5 mg/mL in the presence of a carrier protein (e.g., 0.1% bovine serum albumin) and 0.02-0.05% sodium azide.
Note2: The Ab/ PE-iFluor647 conjugate solution could be stored at 4 °C for two months and kept from light.

Disclaimer: This product is for research use only and is not intended for therapeutic or diagnostic applications. Please contact our technical service representative for more information.