



Rabbit Anti-T7 conjugated to DyLight® 650

Product Number D10-1832

Amount 100 µg total protein

Form/ Storage

Conjugate supplied as a liquid. Upon receipt, store at 2-8°C in the dark.

Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. Avoid exposure to heat and light. Do not freeze.

Buffer

The product is in 100 mM sodium phosphate (pH 7.4), 50 mM sucrose, 0.1% BSA as a stabilizer, and 2 mM sodium azide as a preservative.

Stability

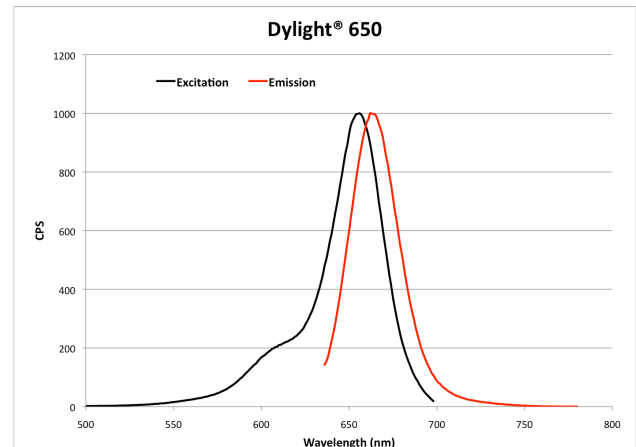
Product should be stored at 2-8°C in the dark and be used within 12 months. If further dilution of the conjugate is required, use diluted material within one week.

Note

For research use only, not for diagnostic or therapeutic use.

Antigen Info

Antibody was raised against T7-tag with amino acid sequence MASMTGGQMG.



Spectral Characteristics

Visible absorption maxima 652

Emission maximum 672

Fluor:Protein ~4

References

Itoh T, Fujita N, Kanno E, Yamamoto A, Yoshimori T, Fukuda M. Golgi-resident Small GTPase Rab33B Interacts with Atg16L and Modulates Autophagosome Formation. *Mol Biol Cell*. 2008 Jul; 19(7): 2916-2925.

Day B, Dahlbeck D, Staskawicz BJ. NDR1 Interaction with RIN4 Mediates the Differential Activation of Multiple Disease Resistance Pathways in Arabidopsis. *Plant Cell*. 2006 Oct; 18(10): 2782-2791.

Rayala SK, Martin E, Sharina IG, Molli PR, Wang X, Jacobson R, Murad F, Kumar R. Dynamic interplay between nitration and phosphorylation of tubulin cofactor B in the control of microtubule dynamics. *PNAS*. 2007 Dec 4; 104(49): 19470-19475.



COLUMBIA BIOSCIENCES