



Rabbit anti-T7 tag IgG conjugated to Dylight ® 650

Product Number D10-1832
Lot Number XXXX
Approximately 5-7
Amount 100 µg
Concentration 1 mg/mL
Store at 2-8°C

Immunogen

T7 (MASMTGGQMG)

Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at -20°C in the dark. Upon Reconstitution store at 2-8°C. Phycobiliproteins are sensitive to freeze-thaw cycles.

Handling

We recommend that the investigator determine the appropriate working concentration for their specific application. Avoid exposure to heat and light.

Buffer

100mM NaPO₄ (pH 7.4), 100mM NaCl, 2mM Sodium Azide

Stability

Upon Reconstitution product should be stored at 2-8 °C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

Antigen Info

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

Note

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

Spectral Characteristics

Visible absorption maxima 652
Emission maximum 672

Fluor:Protein = -2.18

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

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com