



Rabbit anti-T 7 tag conjugated to Dylight® 550

Product Number D9-1832
Amount 0.1 mg total protein
Antigen Amino acid sequence MASMTGGQQMG
Store at 2-8°C

Form/Shipping & Storage

Supplied liquid at 1mg/mL. Upon receipt, store at 2-8 °C. Fluorescent dyes are sensitive to light. Please store in the dark and avoid exposure to sunlight.

Handling

We recommend that the investigator determine the appropriate working concentration for their specific application.

Buffer

0.1M sodium phosphate, 0.15M NaCl, pH7.4, 1% BSA, 0.02% NaN₃

Stability

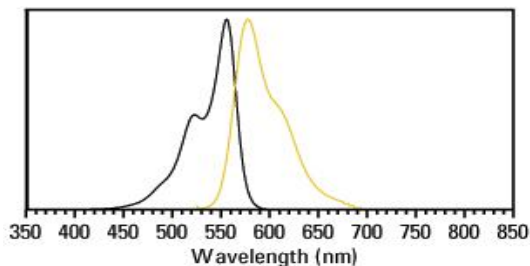
Product should be stored at 2-8°C in the dark and be used within 1 year. If further dilution of the conjugate is required, use diluted material within one month.

Note

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

Spectral Characteristic

Visible absorption maxima 562 nm
Emission maximum 576 nm



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