



## Mouse Anti-6xHistidine (clone AD1.1.10) conjugated to DyLight® 650

Product Number D10-1711-100  
Amount 100 µg  
Store at 2-8°C

### Form/Shipping & 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

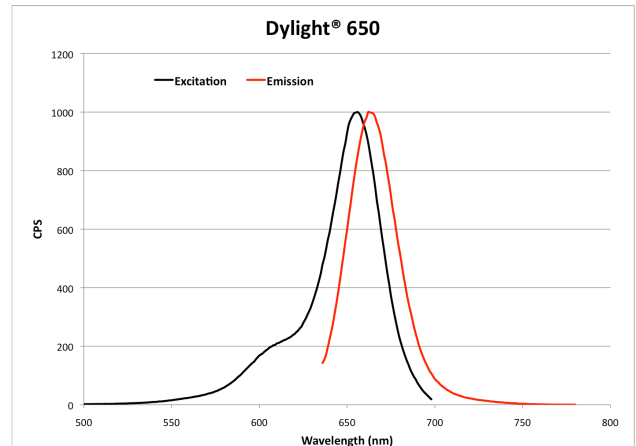
The product is in 10 mM tris (pH 8.2) + 100 mM NaCl + 0.01% NaN<sub>3</sub> + 0.1% BSA.

### 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.



### Spectral Characteristics

Visible absorption maxima 652  
Emission maximum 672

Fluor/IgG ~4

### References

Tang YQ, Han SY, Zheng H, Wu L, Ueda M, Wang XN, Lin Y. Construction of cell surface-engineered yeasts displaying antigen to detect antibodies by immunofluorescence and yeast-ELISA. Appl Microbiol Biotechnol. 2008 Jul;79(6):1019-26.

Kojima T, Takei Y, Ohtsuka M, Kawarasaki Y, Yamane T, Nakano H. PCR amplification from single DNA molecules on magnetic beads in emulsion: application for high-throughput screening of transcription factor targets. Nucleic Acids Res. 2005 Oct 6;33(17):e150.

Tran TM, Moreno A, Yazdani SS, Chitnis CE, Barnwell JW, Galinski MR. Detection of a Plasmodium vivax erythrocyte binding protein by flow cytometry. Cytometry A. 2005;63(1):59-66.