

Mouse Anti-DYKDDDDK IgG conjugated to PerCP

Product Number D11-1718
Amount 100 µg total protein
Store at 2-8°C

Form/Shipping & Storage

Supplied lyophilized. Upon receipt, store at -20 °C. Reconstitute with 1 ml of ddH₂O and store at 2-8 °C. PerCP is 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

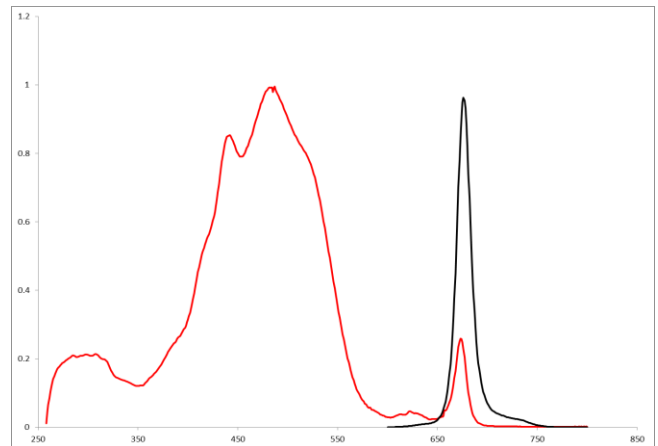
Upon reconstitution, the product is in 10 mM tris (pH 8.2) + 150 mM NaCl + 0.01% NaN₃ + 0.1% BSA. The concentration is 0.1 mg/mL

Stability

Product should be stored at 2-8°C in the dark and be used within 6 months. 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 Characteristics

Visible absorption maxima	483
Emission maximum	673



COLUMBIA BIOSCIENCES