



Rabbit anti-v5 tag conjugated to Dylight® 550

Product Number D9-1834
Lot# XXXX
Approximately 5-7
Amount 0.1 mg total protein
Antigen Amino acid sequence GKPIPPLLGLDST
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, 0.1M NaCl, pH7.4 and 2mM 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.

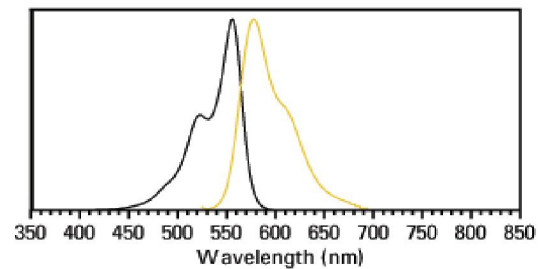
Note

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

Spectral Characteristic

Visible absorption maxima 562 nm

Emission maximum 576 nm



References

Brown GD, Herre J, Williams DL, Willment JA, Marshall ASJ, Gordon S. Dectin-1 Mediates the Biological Effects of β -Glucans. *J. Exp. Med.* 2003 May 5; 197(9):119–1124.

Six E, Ndiaye D, Laabi Y, Brou C, Gupta-Rossi N, Israel A, Logeat F. The Notch ligand Delta1 is sequentially cleaved by an ADAM protease and γ -secretase. *PNAS.* 2003 Jun 24;100(13):7638–7643.