



Rabbit Anti-EP4 Receptor (C-term.) conjugated to SureLight® APC

Product Number D3-1866
Amount 50 µg total protein

Form/ Storage

Conjugate supplied as a lyophilized powder. Upon receipt, store at 2-8°C in the dark.

Handling

Avoid exposure to heat and light. Prior to use reconstitute to 0.5 ml with distilled deionized water, vortex and allow it to sit on ice for 20 minutes.

Buffer

Upon reconstitution, the product is in 10 mM sodium tris (pH 8.2), 50 mM sucrose, 150 mM sodium chloride, 0.1% BSA, 1 µg/mL pentachlorophenol. The concentration of the conjugate is 100 µg/ml.

Stability

Lyophilized material is stable for one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 6 months.

Note

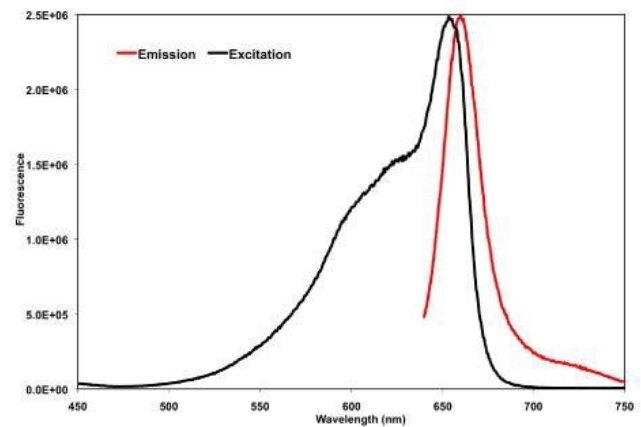
For research use only, not for diagnostic or therapeutic use.

Antigen Info

EP4 receptor C-terminal amino acids 459-488 (GSGRAGPAPKGSLSLQVTFPSETLNLSEKCI)

References

- An, S., Yang, J., Xia, M., et al. Cloning and expression of the EP2 subtype of human receptors for prostaglandin E2. *Biochem. Biophys. Res. Commun.* 197, 263-270 (1993).
- Bastien, L., Sawyer, N., Grygorczyk, R., et al. Cloning, functional expression, and characterization of the human prostaglandin E2 receptor EP2 subtype. *J. Biol. Chem.* 269, 11873-11877 (1994).
- Coleman, R.A., Eglén, R.M., Jones, R.L., et al. Classification of prostanoid receptors IUPHAR receptor compendium. *IUPHAR Compendium* 1-12 (1997).
- Narumiya, S., Sugimoto, Y., and Ushikubi, F. Prostanoid receptors: structures, properties, and functions. *Physiol. Rev.* 79, 1193-1226 (1999).



Fluorescence excitation and emission spectra of APC in 100 mM sodium phosphate (pH 7.2) + 1 mM EDTA and 1 mM sodium azide. The emission scan was taken with excitation at 630 nm. The excitation scan was taken with an emission at 660 nm. The curves were normalized to equalize peak heights.

Spectral Characteristics

Visible absorption maxima 650
Emission maximum 657 nm

Concentration upon Reconstitution

SureLight APC= 0.49 µM
α-EP4= 0.33 µM

Reactivity

Human, murine, rat, and ovine EP4 receptor; non-reactive with EP1, EP2, and EP3 receptors; other species not tested.

4985 Winchester Blvd. Frederick, MD 21703

301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com

For sales inquiries: sales@columbiabiosciences.com