



## Goat anti-Glutathione-S-Transferase IgG conjugated to SureLight® P-3

Product Number D7-1310  
Amount 100 µg  
Store at 4°C

### Form/Shipping & Storage

Supplied as a lyophilized powder. Upon receipt, store at 4°C in the dark. Phycobiliproteins are sensitive to freeze-thaw cycles; after reconstitution, store at 2-8°C in the dark, do not freeze.

### Handling

Prior to use, reconstitute to 200 µl with distilled deionized water, vortex gently and allow vial to sit on ice for 20 minutes. Avoid exposure to heat and light.

### Buffer

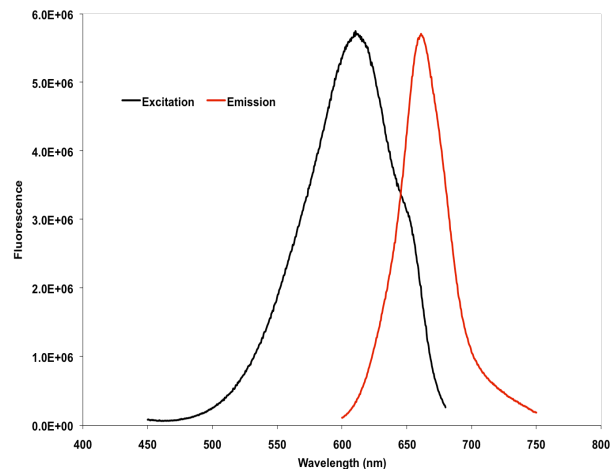
Upon rehydration with 200 µl distilled deionized water, the product is in 100 mM sodium phosphate (pH 7.4), 100 mM sucrose, 150 mM sodium chloride, 0.1% BSA as a stabilizer, and 2 mM sodium azide as a preservative.

### Stability

Lyophilized material is stable for up to one year. After product has been reconstituted, product should be stored at 2-8°C in the dark and be used within 3 months. If further dilution of the conjugate is required, use diluted material within one week.

### Use Restriction

Buyer hereby agrees to the following use restriction: Buyer will not use or sell the SureLight® P1 or P3 detection reagent purchased from Columbia Biosciences for the production, use or sale of multi-well plates or multi-chamber slides containing at least four (4) wells or chambers per test device, with each well or chamber containing an array of immobilized proteins, peptides or nucleic acids in research applications of multi-well arrays that do not require approval by state or federal regulatory agencies. Such rights are currently licensed and are the sole property of the licensee.



### Note

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

### Spectral Characteristics

Visible absorption maxima	614>650 (sh) nm
Emission maximum	662 nm



**COLUMBIA BIOSCIENCES**